

Tired of ELISA? – It's Time for Lumit™ Immunoassays

An Easier and Faster Method for Protein Detection

Welcome

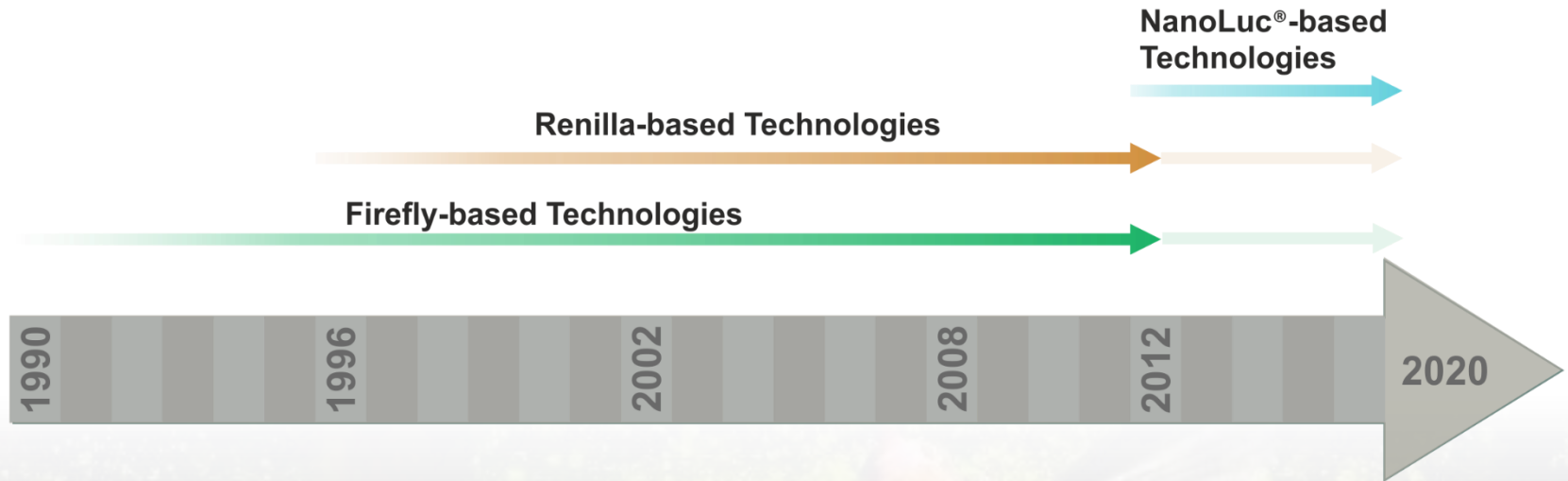
Dr. Erik Bonke | Application Specialist | Promega GmbH





Promega – The Bioluminescent Company

A Continuously Grown Expertise in Luciferase-based Technologies



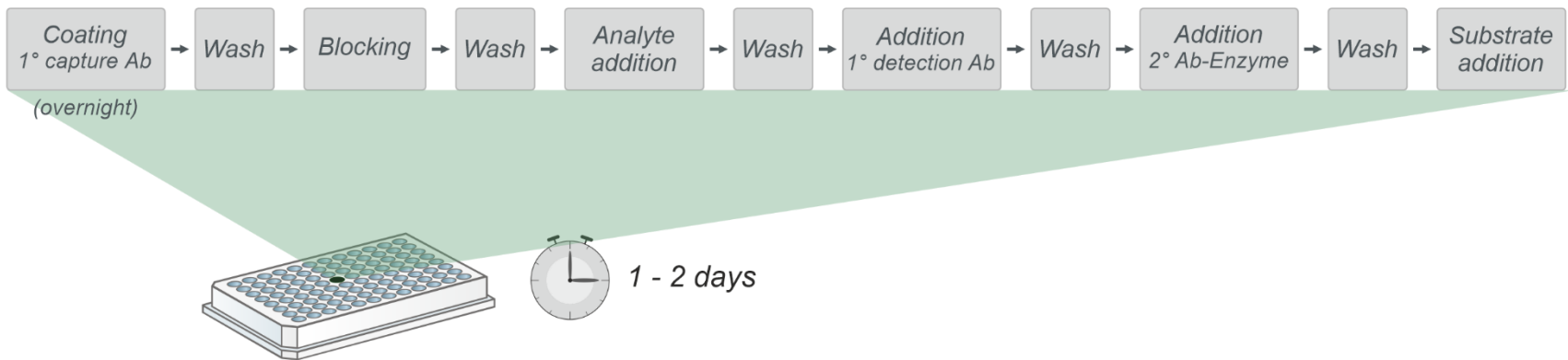
- Reporter Gene Assays
- GloSensor™ (cAMP, Protease Assays)
- GloResponse™ (Signaling Pathways)
- Rapid Response™ (Signaling Pathways)
- Cell-Health Assays
- Bioassays (ADCC, PDL1)
- NanoBRET™ / NanoBiT® (Protein Interaction)
- NanoBRET™ Target Engagement
- HiBiT Protein Tagging System
- Lumit™ Immunoassays



Lumit™ Immunoassays

The Powerful Alternative to Conventional Immunoassay Approaches

Traditional ELISA Workflow



- Traditional ELISA is a heterogenous multistep process involving several wash / incubation steps

Lumit™ Immunoassays

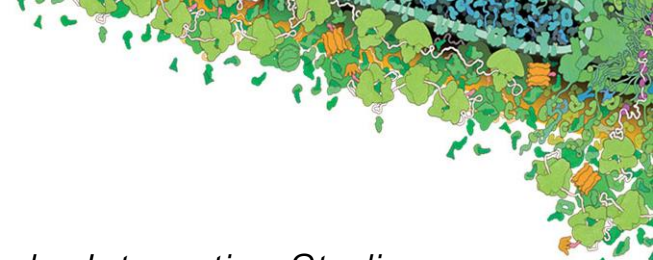
The Powerful Alternative to Conventional Immunoassay Approaches

Traditional ELISA Workflow



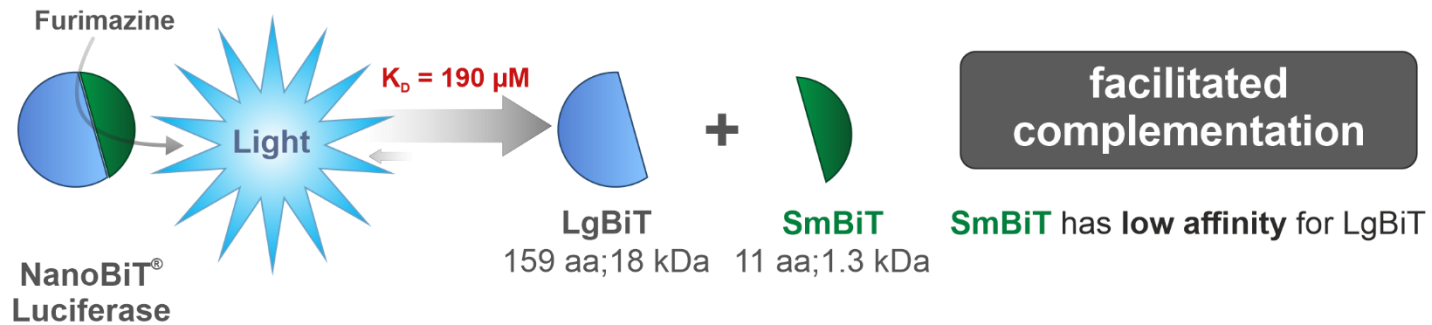
- Traditional ELISA is a heterogenous multistep process involving several wash / incubation steps
- Based on NanoLuc® luciferase we developed *Lumit™ Immunoassays*
 - ✓ Easy and fast (30 – 120 min)
 - ✓ High Sensitivity (low number of cells)
 - ✓ Broad dynamic range (3 – 4 logs)
 - ✓ Flexible formats (96- or 384-well)
 - ✓ Homogenous and HTS compatible

Hwang, B. *et al.* (2020) *Commun Biol.* 3:8



NanoLuc[®] Binary Technology (NanoBiT[®])

A Structural Complementation Reporter Designed for Biomolecular Interaction Studies



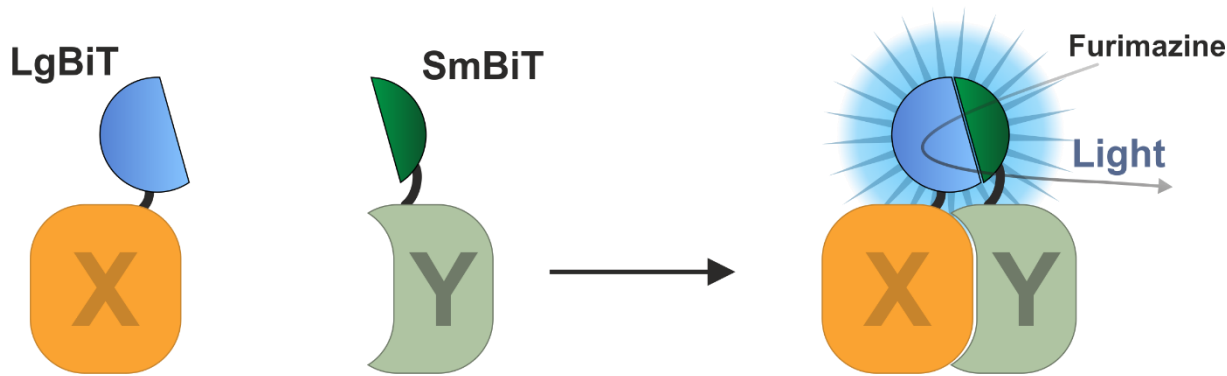
Complementation facilitated through ...

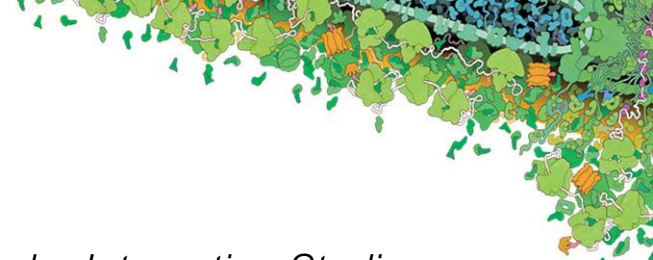
(direct) protein:protein interaction



NanoBiT[®] PPI System

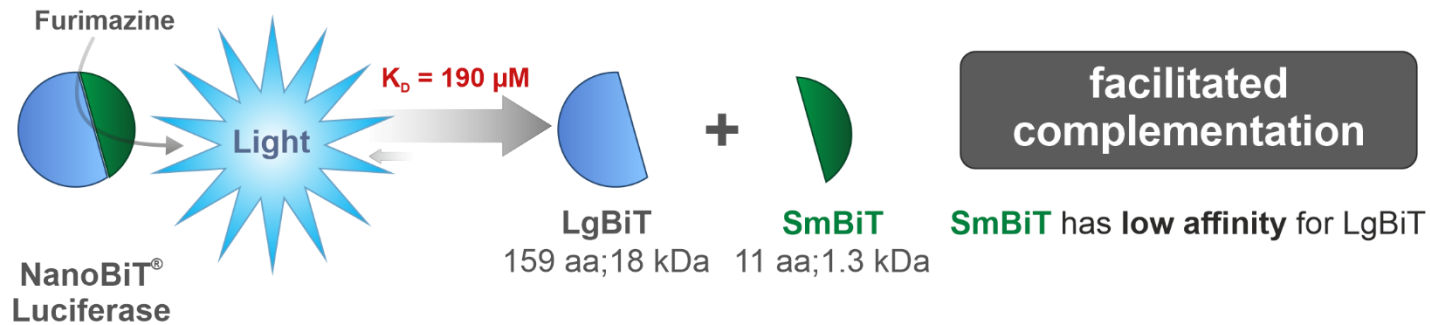
Dixon et al. 2016, ACS Chemical Biology





NanoLuc[®] Binary Technology (NanoBiT[®])

A Structural Complementation Reporter Designed for Biomolecular Interaction Studies



Complementation facilitated through ...

(indirect) Ab:Ab “interaction”



Lumit[™] Immunoassays

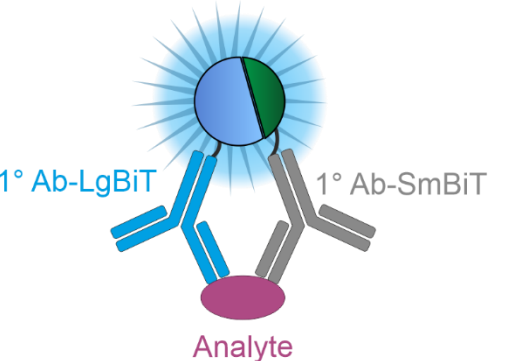
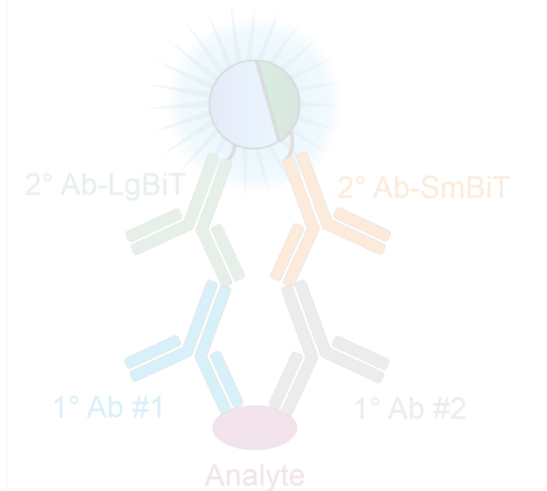
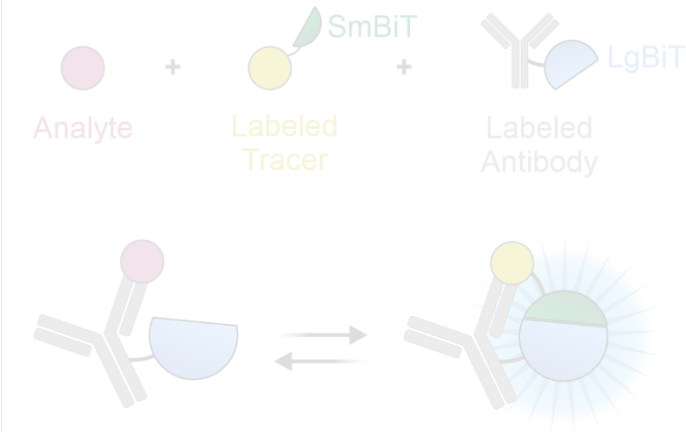
Hwang et al. 2020, Commun Biol.





Lumit™ Immunoassays

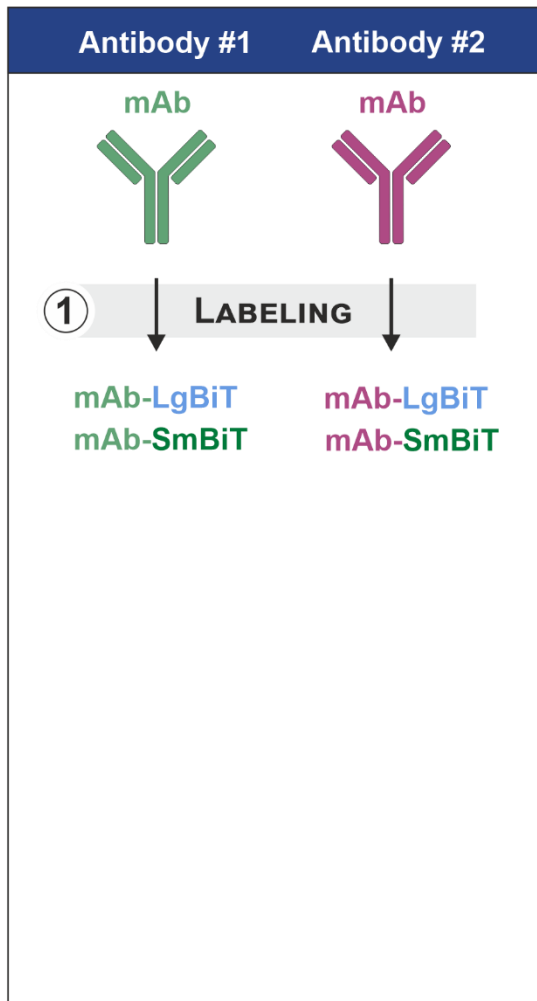
Different Formats for Maximum Flexibility

Direct	Indirect	Competitive
 <p>1° Ab-LgBiT</p> <p>1° Ab-SmBiT</p> <p>Analyte</p>	 <p>2° Ab-LgBiT</p> <p>2° Ab-SmBiT</p> <p>1° Ab #1</p> <p>1° Ab #2</p> <p>Analyte</p>	 <p>Analyte</p> <p>Labeled Tracer</p> <p>Labeled Antibody</p>
<ul style="list-style-type: none">• Requires labeling of 1°Abs• Validated for cytokines and peptide hormones• <i>Ready-to-use</i> assays for<ul style="list-style-type: none">✓ IL1-β, IFN-γ, IL-2, IL-6, IL-10, IL-4, TNF-α, VEGF, ...✓ Insulin and glucagon	<ul style="list-style-type: none">• Avoids labeling of 1°Abs• Generic pre-labeled 2°Abs (different species available)• Validated for intracellular PTMs, e.g. phosphorylation	<ul style="list-style-type: none">• Requires antibody and tracer labeling• Establish competitive antibody binding assays• <i>Ready-to-use</i> assay for hFcRn:Fc protein interactions, e.g. antibodies<ul style="list-style-type: none">✓ Lumit™ FcRn Binding Immunoassay

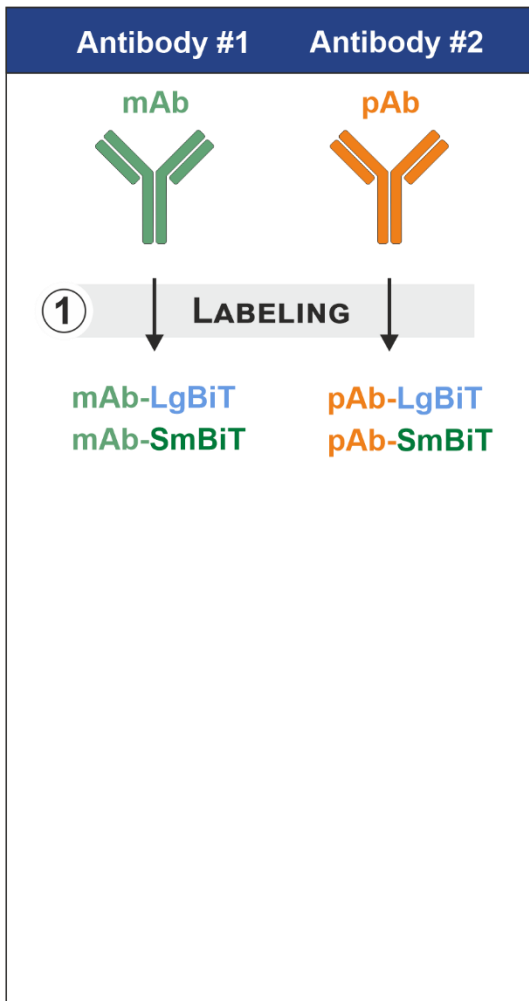


Build-Your-Own Direct Lumit™ Immunoassay

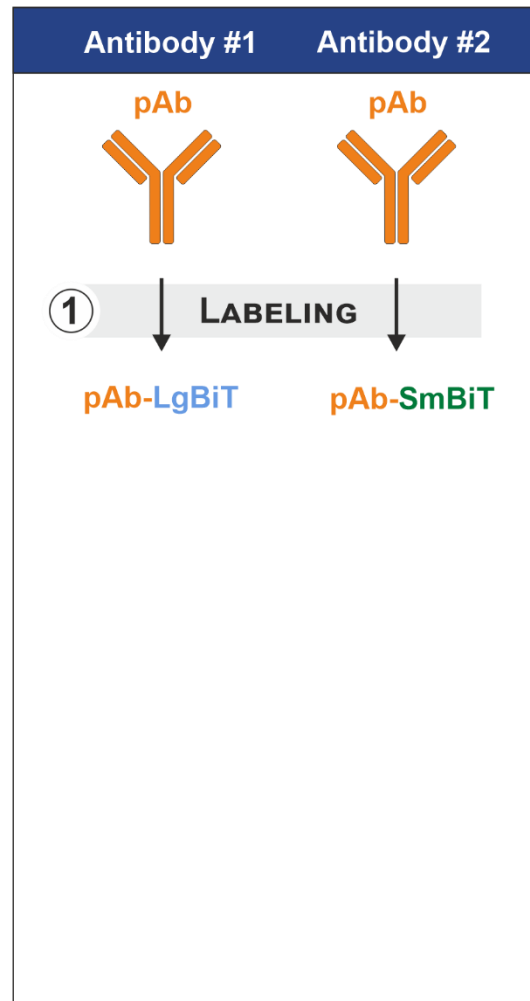
Various Options for Your Convenience



😊 highest specificity and sensitivity



😊 only one mAb needed

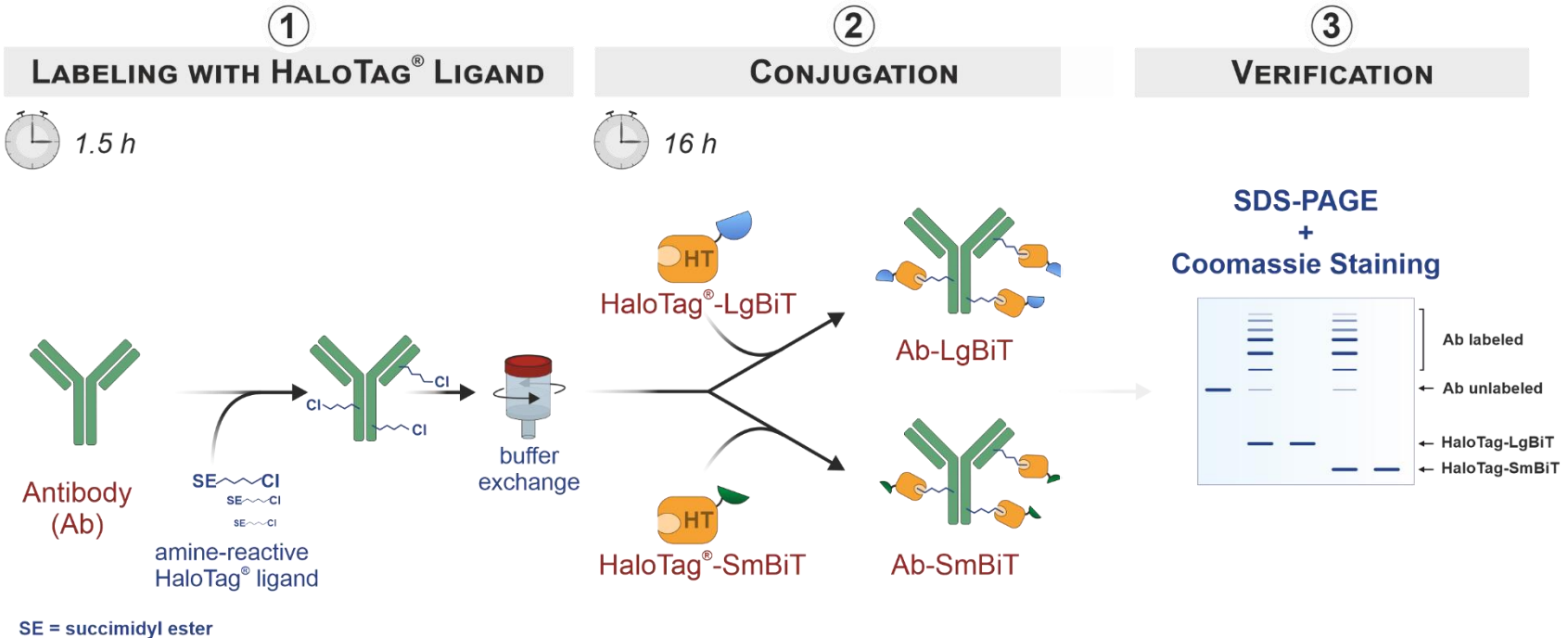


😊 most inexpensive variant



Build-Your-Own Direct Lumit™ Immunoassay

Step 1: Labeling of Antibodies with the Lumit™ Immunoassay Labeling Kit



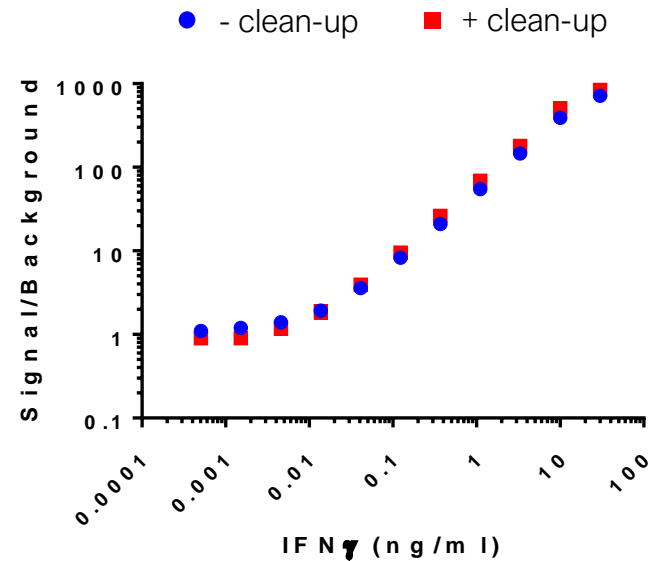
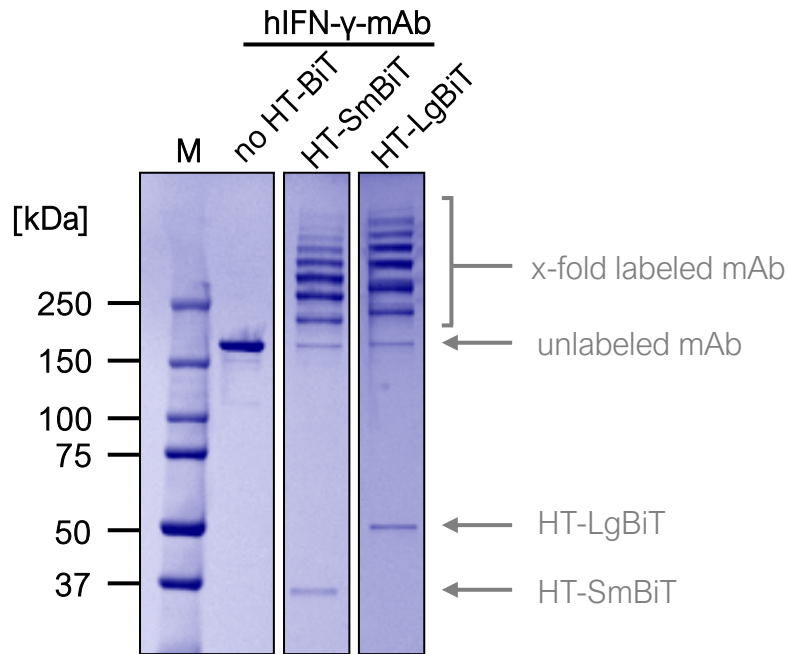
FACTS

- Easy and robust 2-day protocol
- Attachment is highly efficient (> 90%)
- Oriented BiT subunits for maximum activity



Build-Your-Own Direct Lumit™ Immunoassay

Step 1: Labeling of Antibodies with the Lumit™ Immunoassay Labeling Kit



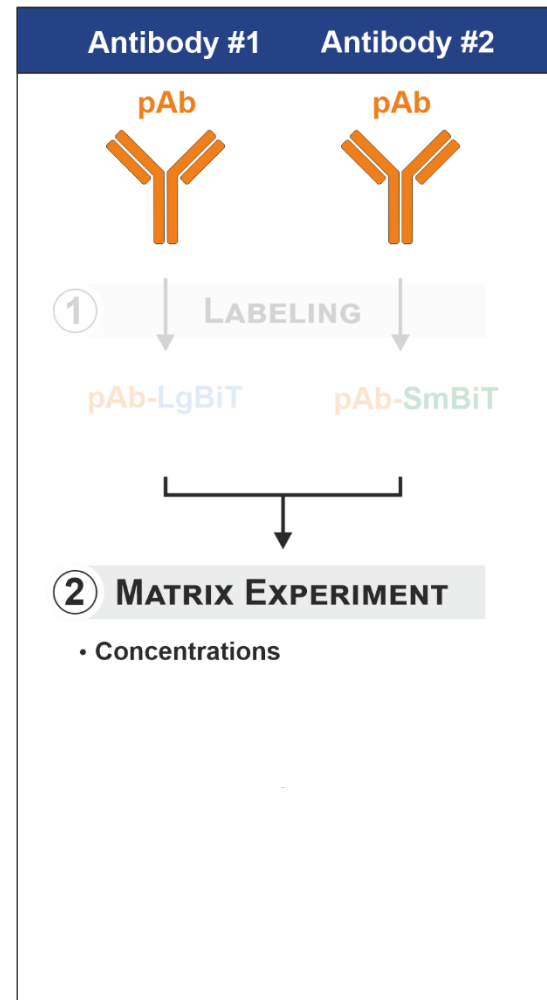
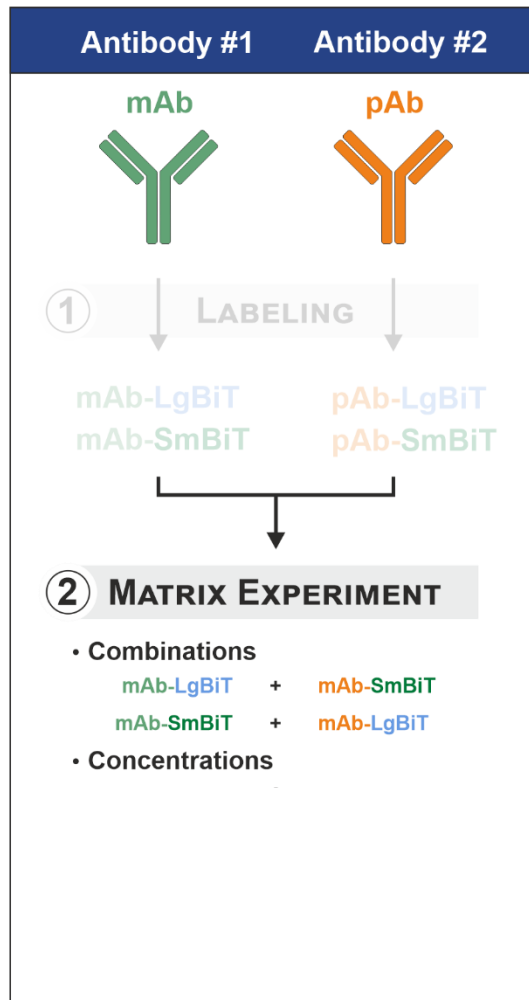
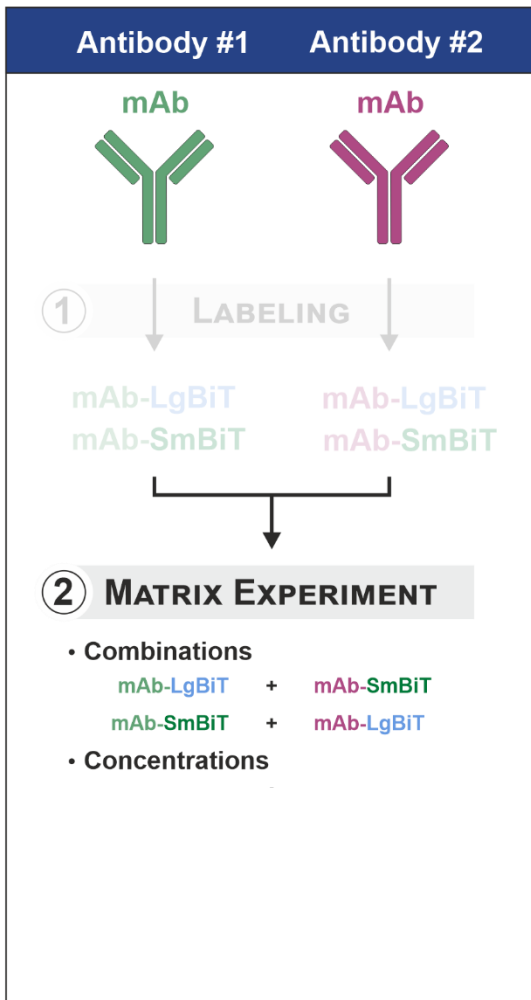
FACTS

- Easy and robust 2-day protocol
- Attachment is highly efficient (> 90%)
- Oriented BiT subunits for maximum activity



Build-Your-Own Direct Lumit™ Immunoassay

Various Options for Your Convenience



😊 highest specificity and sensitivity

😊 only one mAb needed

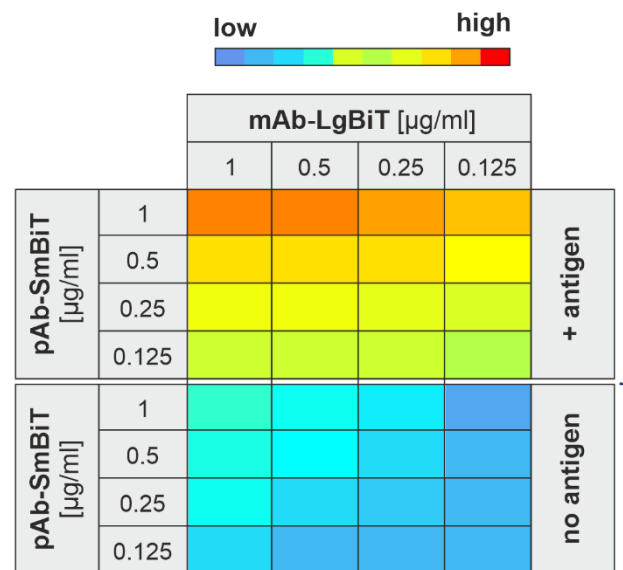
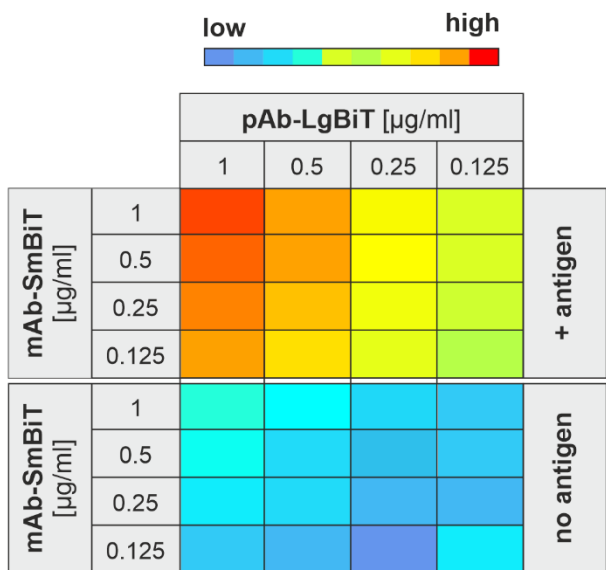
😊 most inexpensive variant



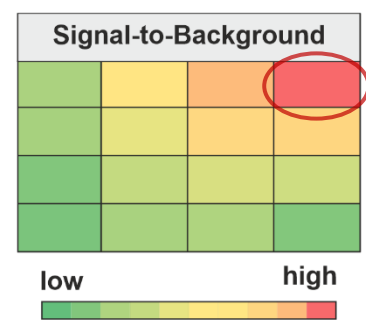
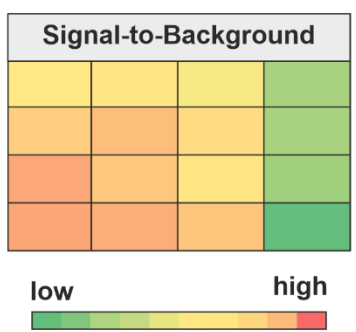
Build-Your-Own Direct Lumit™ Immunoassay

Step 2: Identification of best antibody combination/concentration

Matrix experiment to determine maximal signal-to-background (S/B) ratio



Background

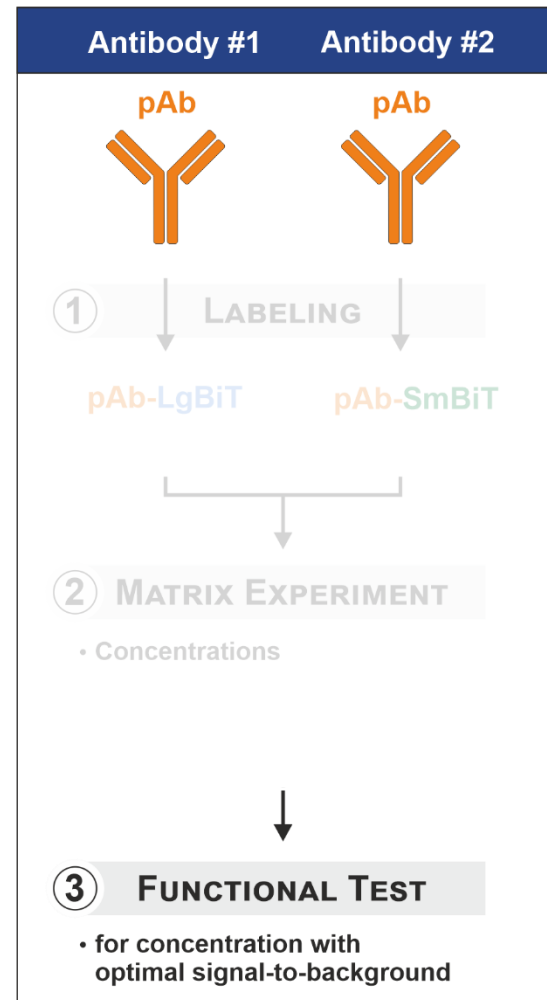
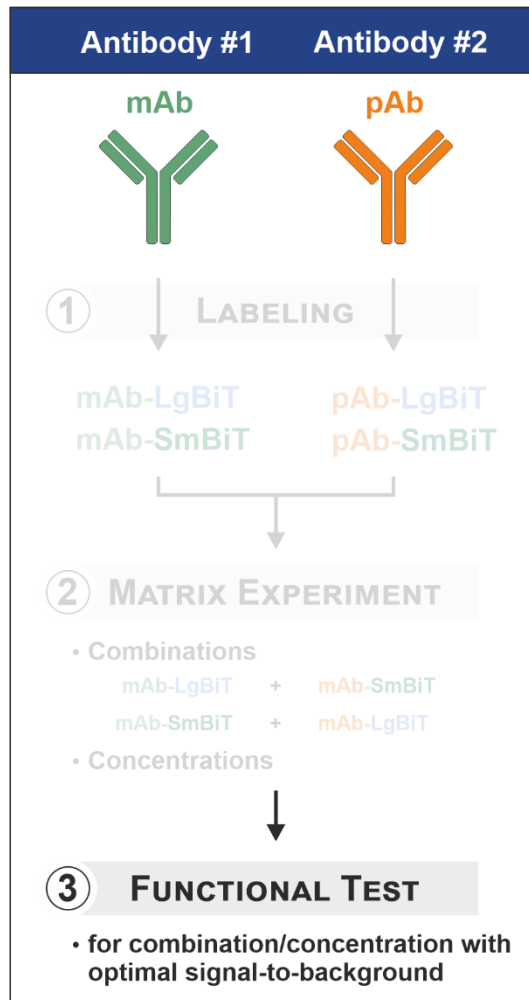
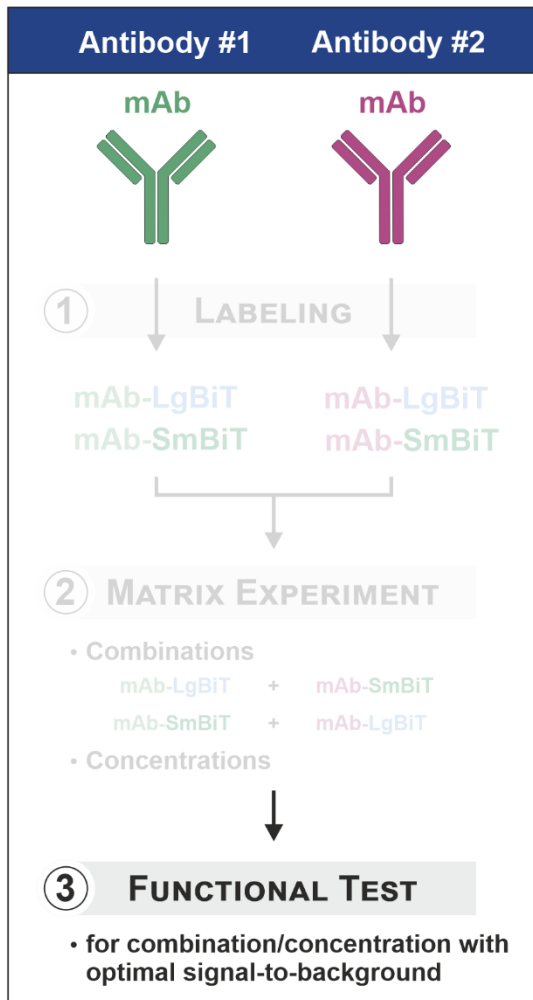


mAb-LgBiT
0.125 $\mu\text{g/ml}$
+
pAb-SmBiT
1 $\mu\text{g/ml}$



Build-Your-Own Direct LumitTM Immunoassay

Various Options for Your Convenience



😊 highest specificity and sensitivity

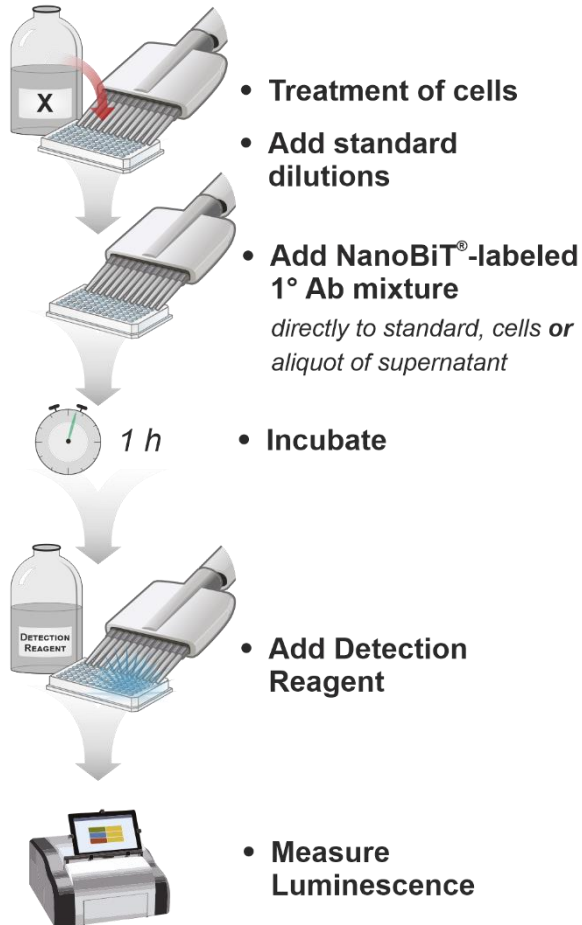
😊 only one mAb needed

😊 most inexpensive variant

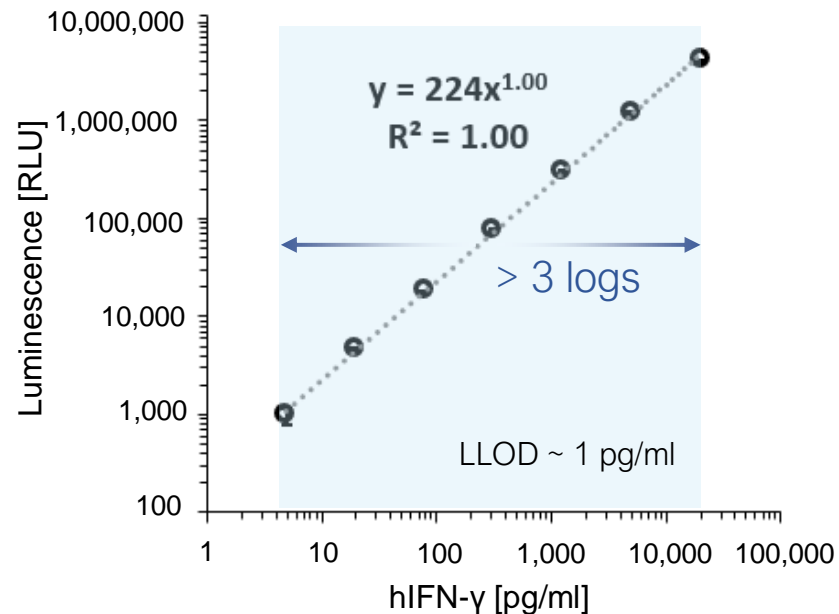


Build-Your-Own Direct Lumit™ Immunoassay

Step 3: Functional Test of Identified Antibody Combination



Lumit™ (human) IFN- γ Immunoassay



OBSERVATIONS

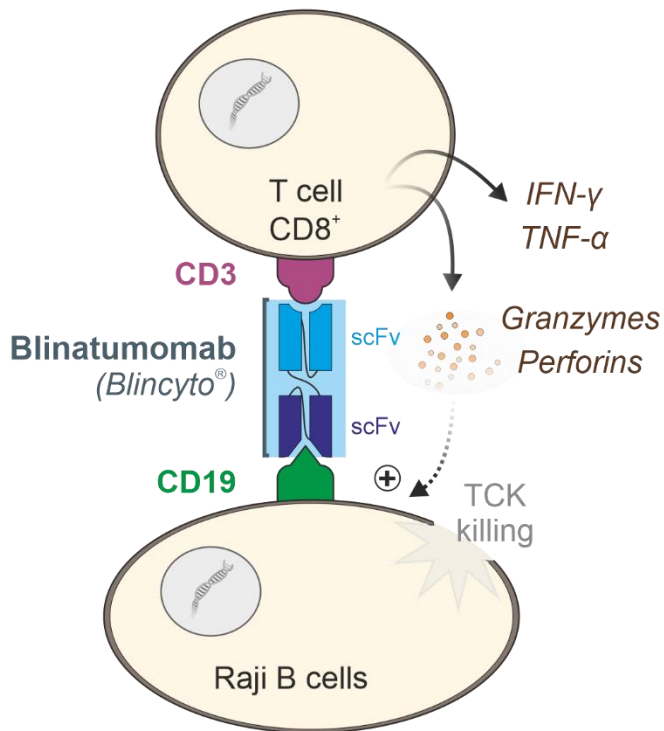
- Good sensitivity and broad dynamic range (≥ 3 logs)
- No need for sample dilution



Build-Your-Own Direct Lumit™ Immunoassay

Step 3: Functional Test of Identified Antibody Combination

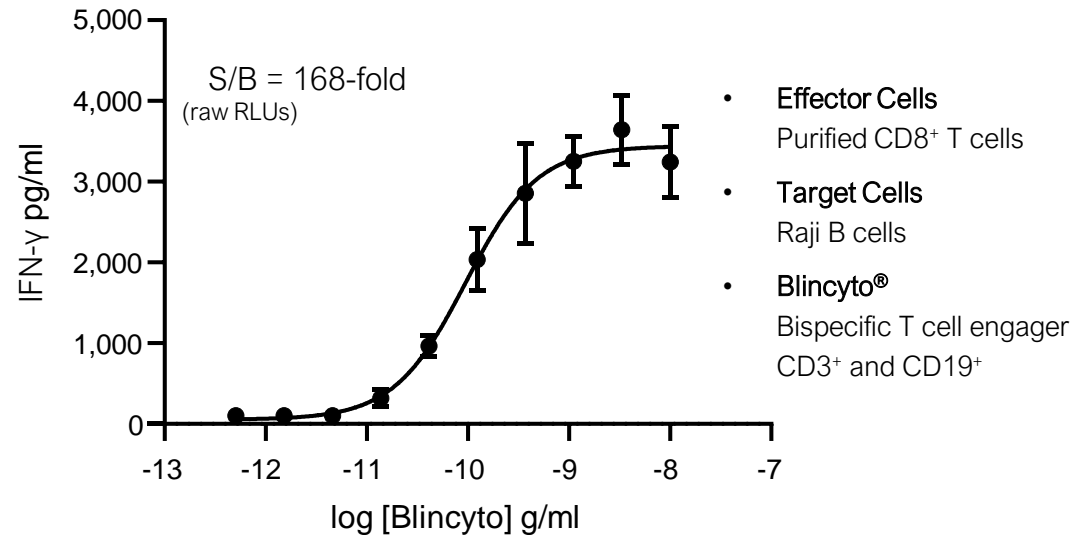
BiTE-induced IFN- γ release from CD8⁺ T cells



BiTE: Bispecific T cell engager

scFv: single-chain variable fragment

Lumit™ (human) IFN- γ Immunoassay



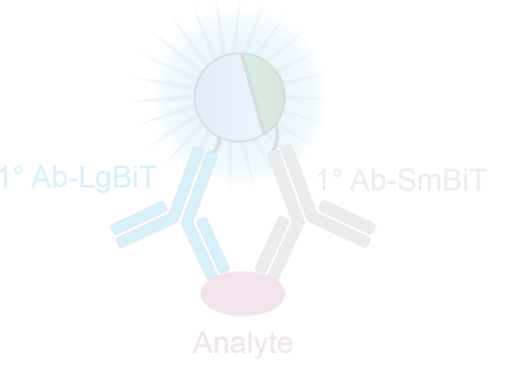
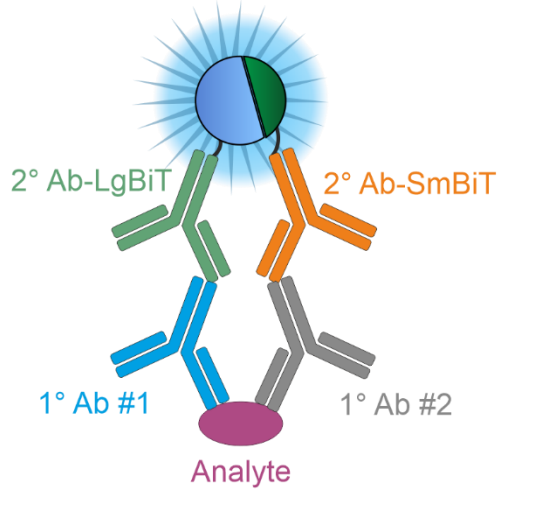
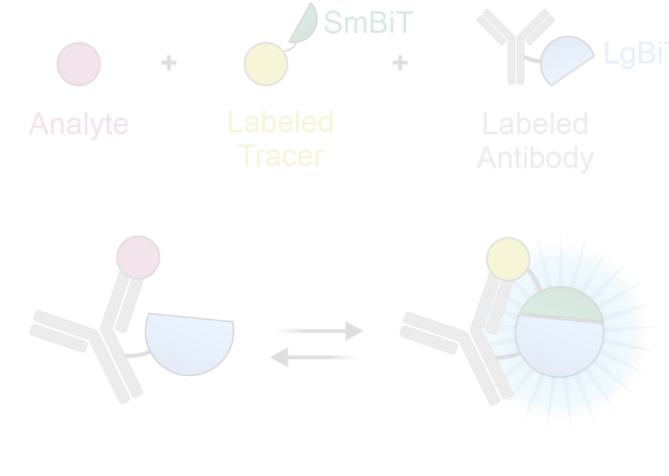
OBSERVATIONS

- Good sensitivity and broad dynamic range (≥ 3 logs)
- No need for sample dilution
- Excellent cell-based performance (addition to cells)



Lumit™ Immunoassays

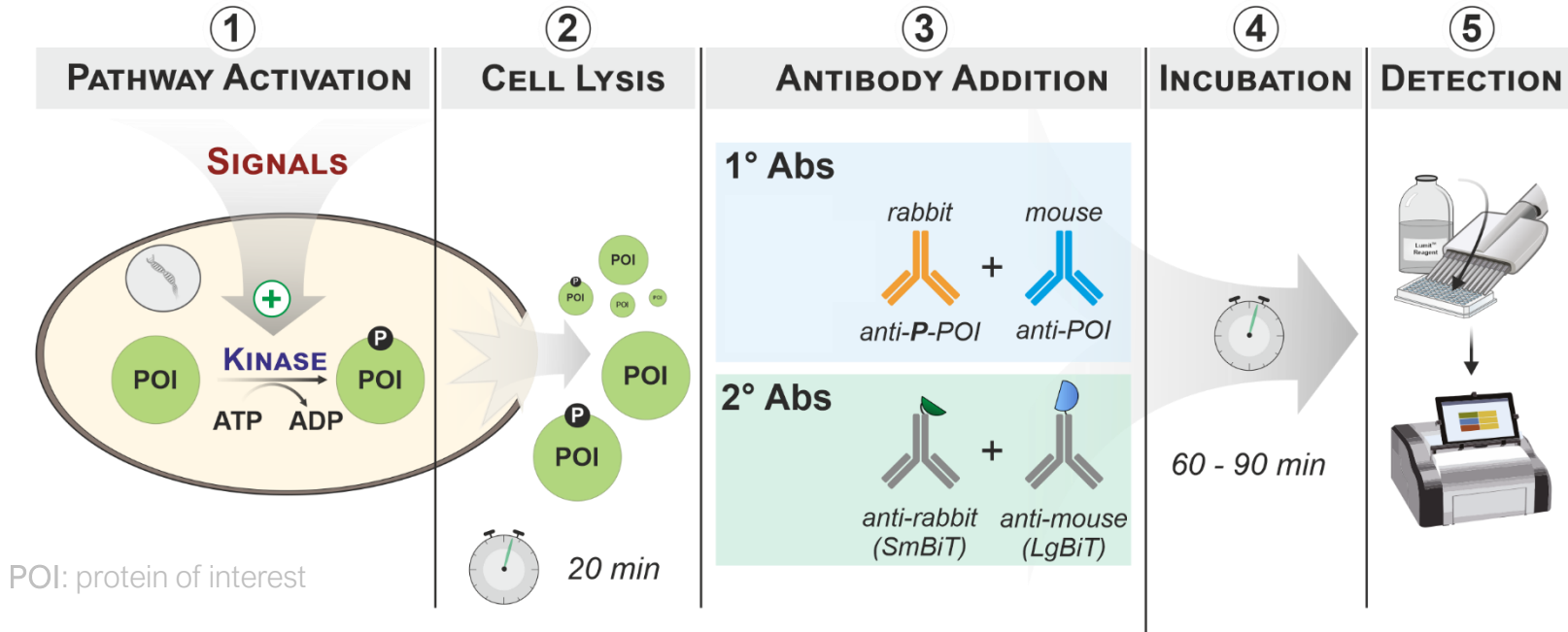
Different Formats for Maximum Flexibility

Direct	Indirect	Competitive
 <p>1° Ab-LgBiT 1° Ab-SmBiT</p> <p>Analyte</p>	 <p>2° Ab-LgBiT 2° Ab-SmBiT</p> <p>1° Ab #1 1° Ab #2</p> <p>Analyte</p>	 <p>Analyte Labeled Tracer Labeled Antibody</p>
<ul style="list-style-type: none">• Requires labeling of 1°Abs• Validated for cytokines and peptide hormones• <i>Ready-to-use</i> assays for<ul style="list-style-type: none">✓ IL1-β, IFN-γ, IL-2, IL-6, IL-10, IL-4, TNF-α, VEGF, ...✓ Insulin and glucagon	<ul style="list-style-type: none">• Avoids labeling of 1°Abs• Generic pre-labeled 2°Abs (different species available)• Validated for intracellular PTMs, e.g. phosphorylation	<ul style="list-style-type: none">• Requires antibody and tracer labeling• Establish competitive antibody binding assays• <i>Ready-to-use</i> assay for hFcRn:Fc protein interactions, e.g. antibodies<ul style="list-style-type: none">✓ Lumit™ FcRn Binding Immunoassay



Lumit™ Immunoassay Cellular Systems

Study Cellular Signaling Events



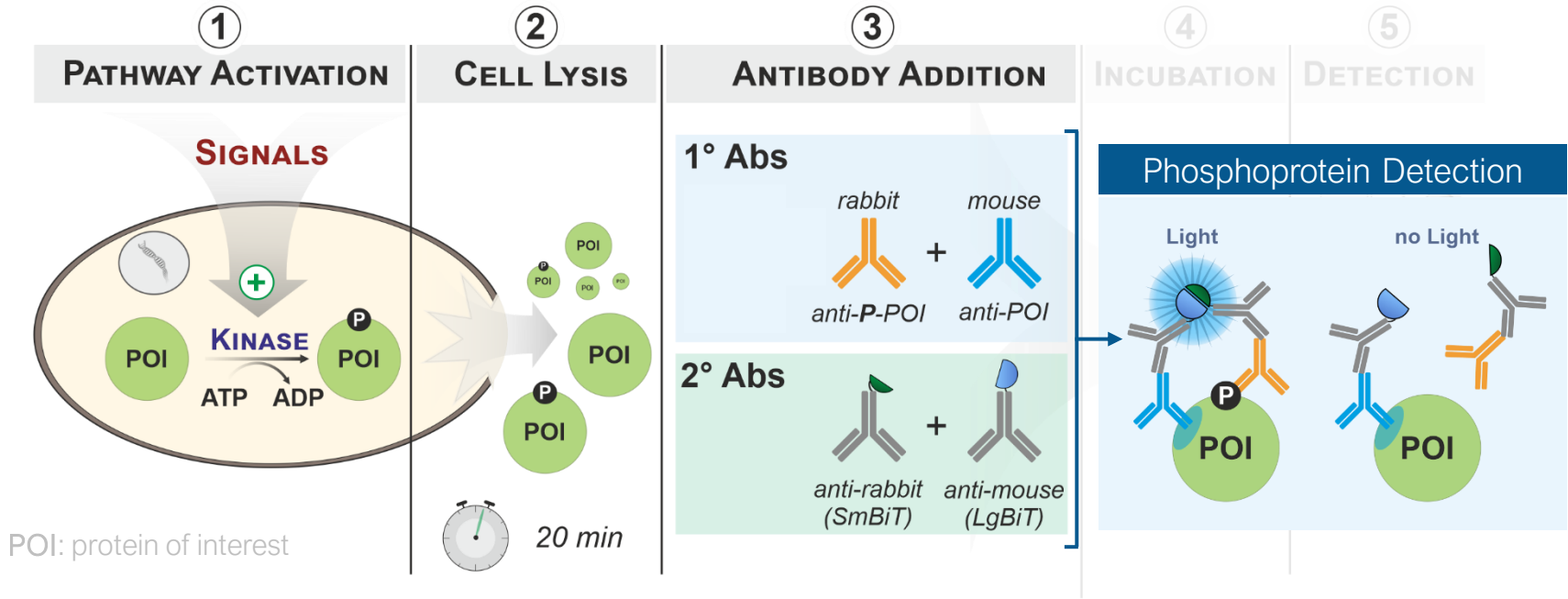
Additionally available pre-labeled 2° Abs:

- anti-rabbit (LgBiT)
- anti-mouse (SmBiT)
- anti-goat (LgBiT)
- anti-goat (SmBiT)



Lumit™ Immunoassay Cellular Systems

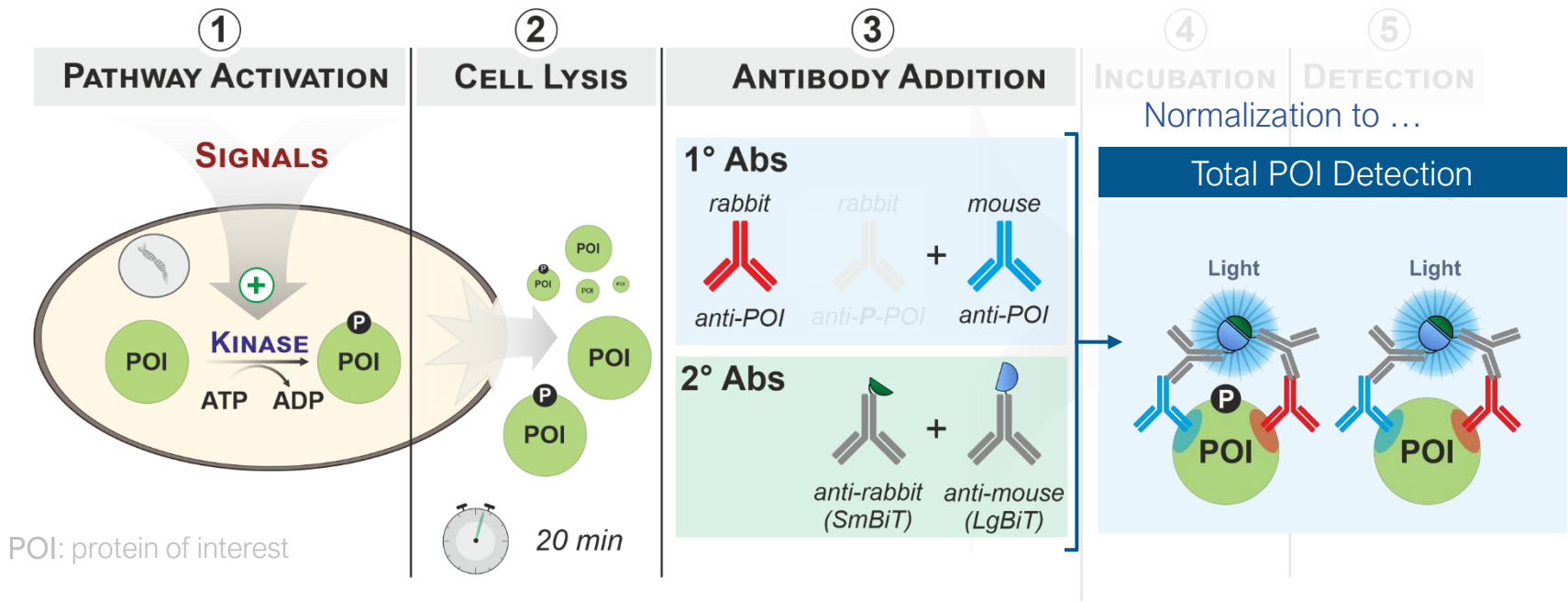
Study Cellular Signaling Events



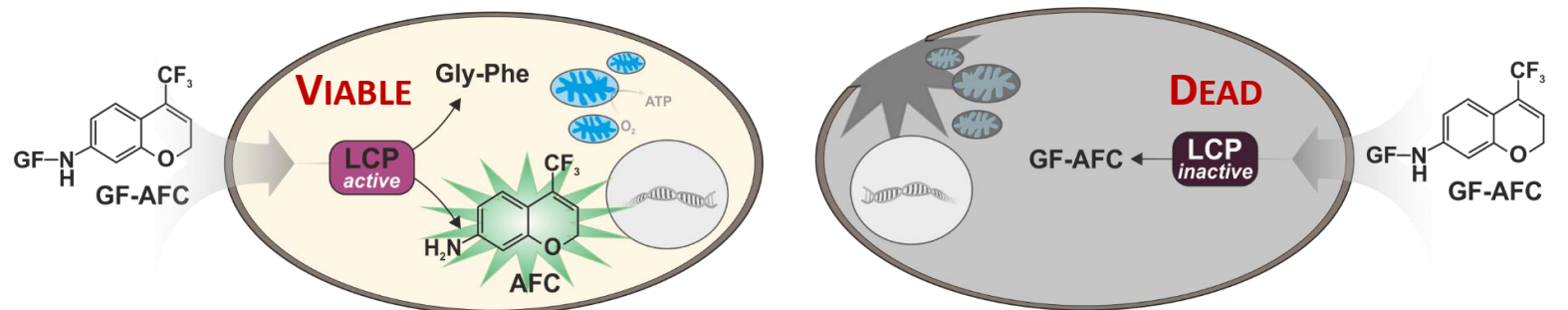


Lumit™ Immunoassay Cellular Systems

Study Cellular Signaling Events



Normalization on number of viable cells using CellTiter-Fluor™

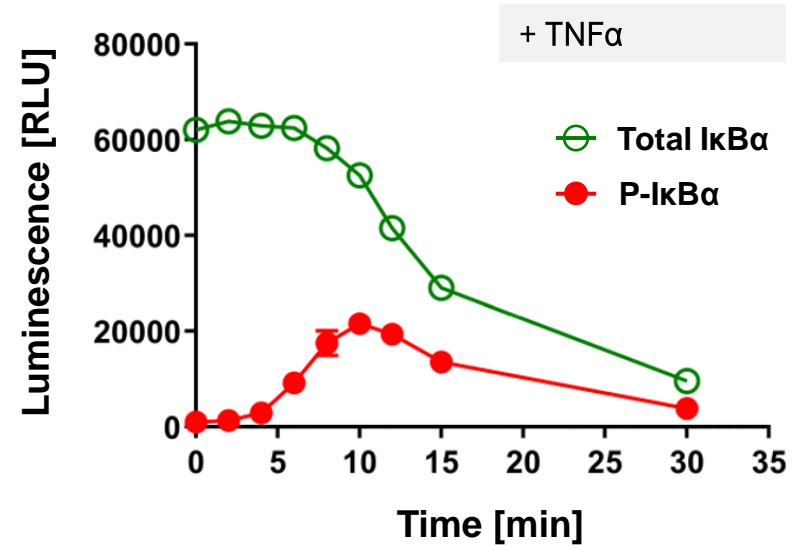
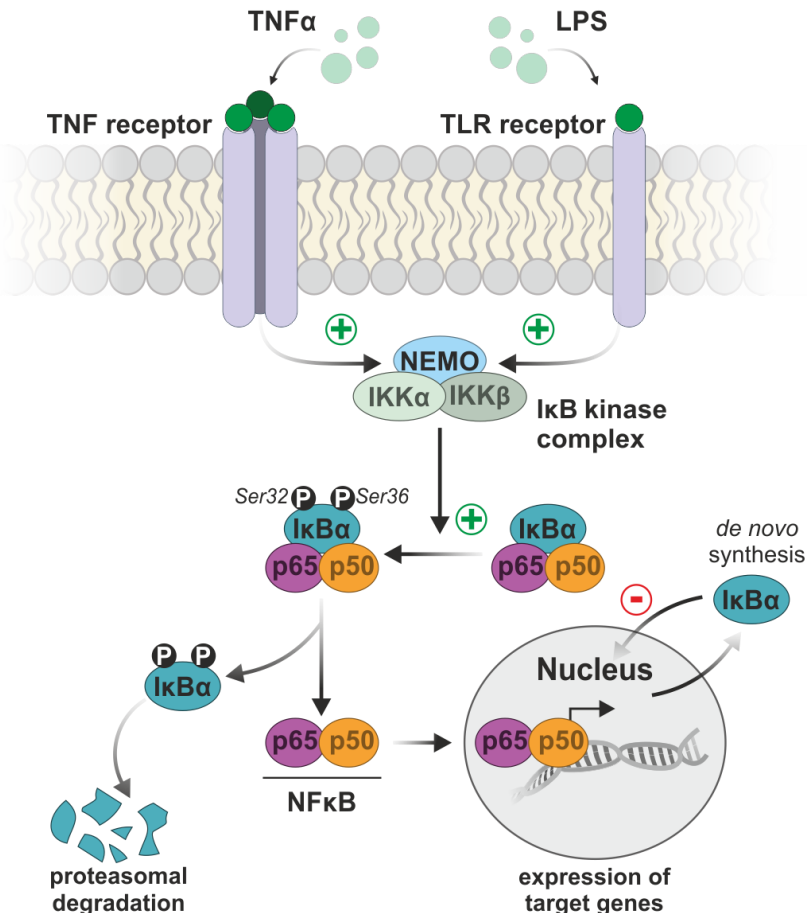


LCP: Live-cell protease



Signaling Pathway and Kinase Activity Analysis

Studying the NFκB Pathway



Modified from Hwang, B. *et al.* (2020) *Commun Biol.* 3:8

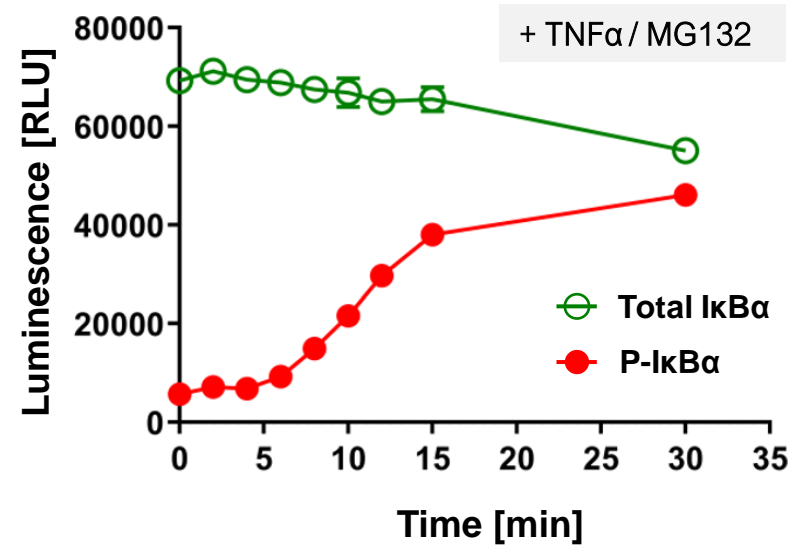
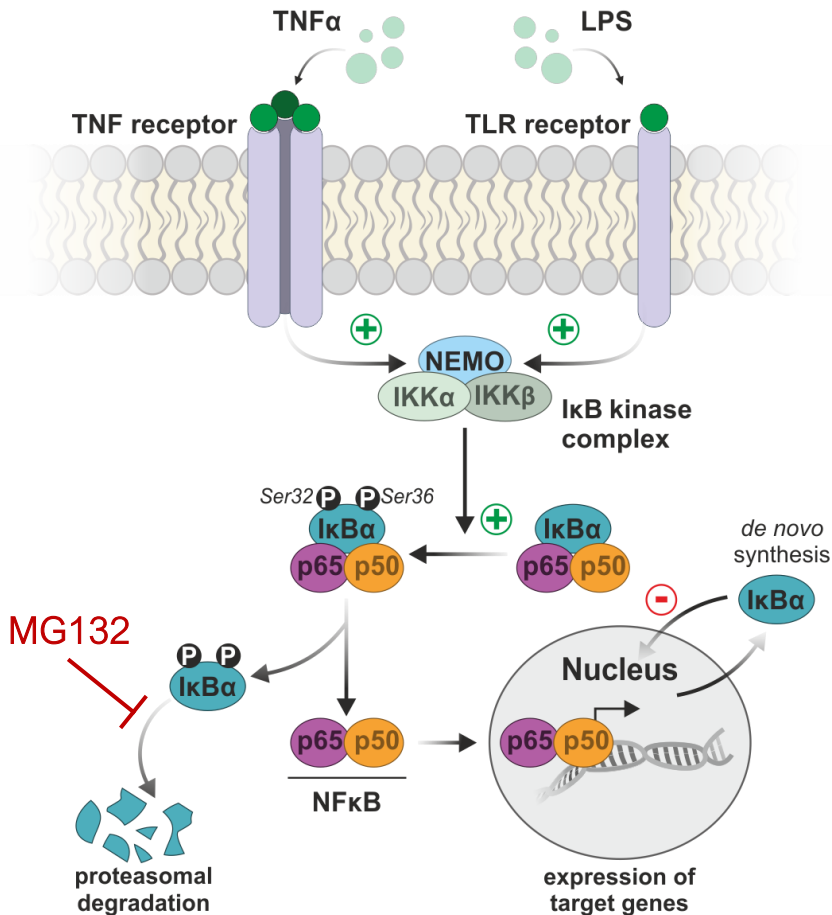
OBSERVATION

- IκBα phosphorylation at Ser32 (pS32)
- Immediately followed by rapid degradation



Signaling Pathway and Kinase Activity Analysis

Studying the NFκB Pathway



Modified from Hwang, B. *et al.* (2020) *Commun Biol.* 3:8

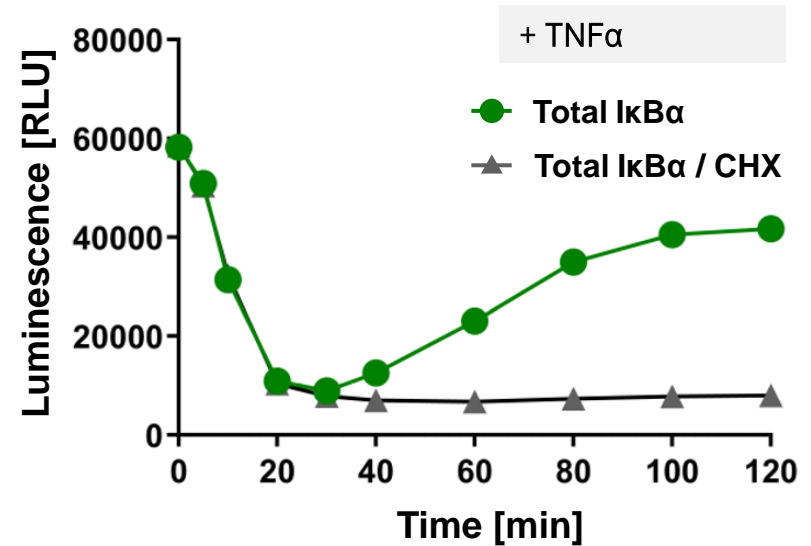
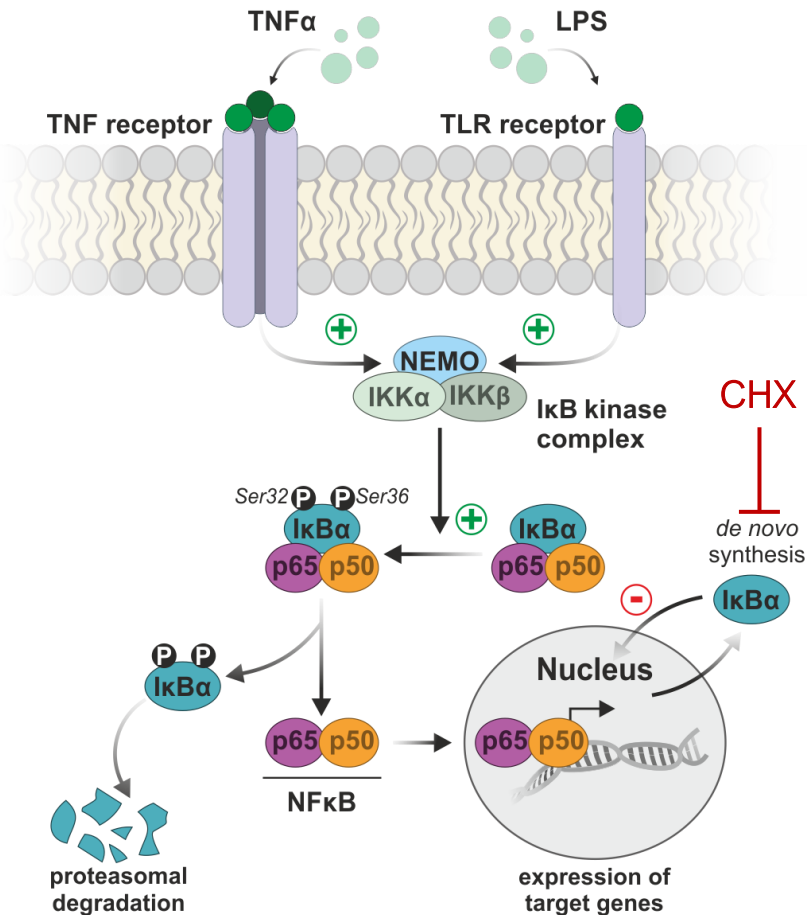
OBSERVATION

- Decrease in IκBα degradation
- Accumulation of phosphorylated IκBα



Signaling Pathway and Kinase Activity Analysis

Studying the NFκB Pathway



Modified from Hwang, B. *et al.* (2020) *Commun Biol.* 3:8

OBSERVATION

- Cycloheximide (CHX) blocks NFκB-triggered *de novo* IκBα biosynthesis

Lumit™ Immunoassay Cellular Systems

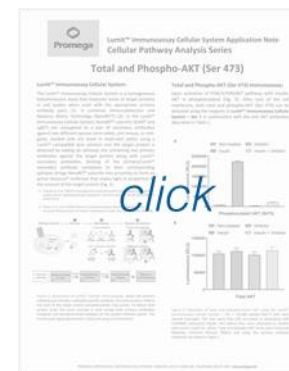
A Universal Immunoassay to Study Cellular Signaling



Validated with >20 phospho- and total proteins using 8 cell types, suggesting this universal immunoassay can be adapted for any pathway **with the appropriate antibodies**

- IκBα (phospho-Ser32 and total protein)
- STAT3 (phospho-Tyr705 and total protein)
- BTK (phospho-Tyr223)
- Estrogen receptor (total protein)
- β-Catenin (total protein)
- CREB (phospho-Ser133)
- P38 MAPK (phospho-Thr180/182)
- NFκB p65 (phospho-Ser536 and total protein)
- AKT (phospho-Ser473, phospho-Thr308, and total protein)
- Retinoblastoma tumor suppressor protein (phospho-Ser807/811 and phospho-Ser780)
- S6 ribosomal protein (phospho-Ser235/236, phospho-Ser240/244)
- MEK1/2 (phospho-Ser217/221, phospho-Ser298)

Application Notes

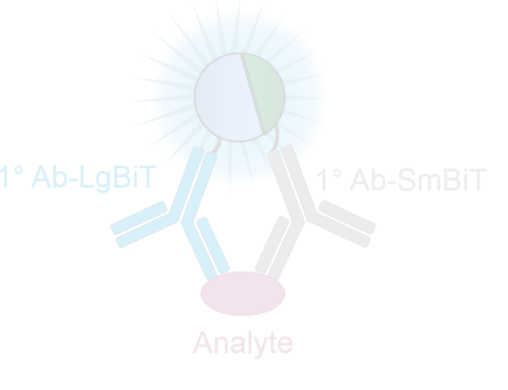
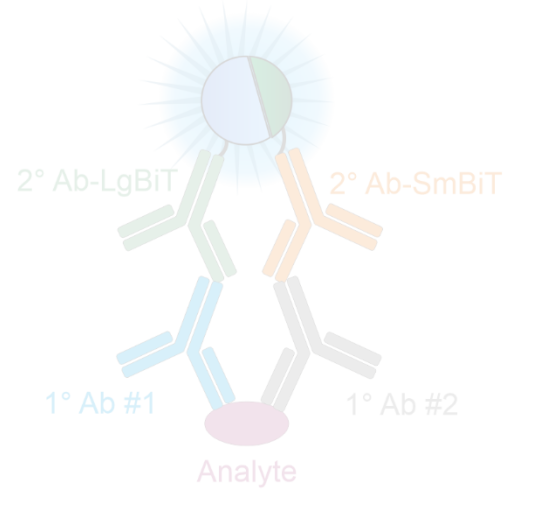
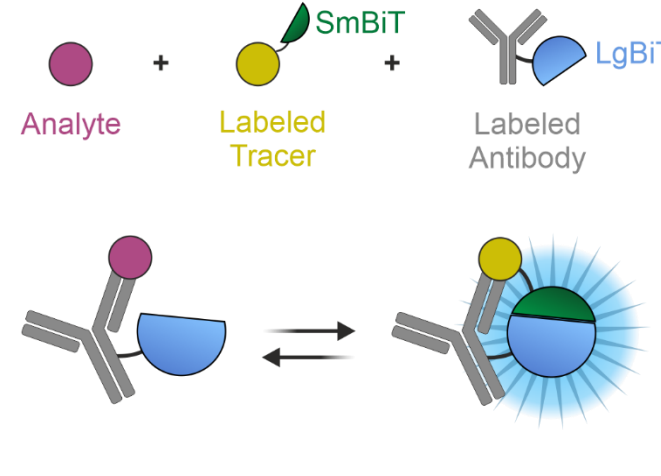


Available on Promega website with information on antibodies used



Lumit™ Immunoassays

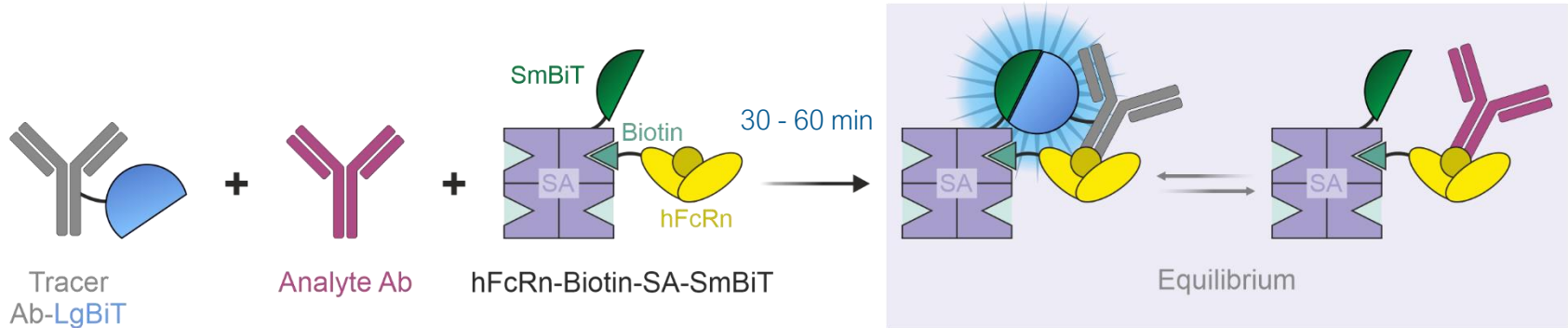
Different Formats for Maximum Flexibility

Direct	Indirect	Competitive
 <p>1° Ab-LgBiT 1° Ab-SmBiT</p> <p>Analyte</p>	 <p>2° Ab-LgBiT 2° Ab-SmBiT</p> <p>1° Ab #1 1° Ab #2</p> <p>Analyte</p>	 <p>Analyte Labeled Tracer Labeled Antibody</p>
<ul style="list-style-type: none"> • Requires labeling of 1°Abs • Validated for cytokines and peptide hormones • <i>Ready-to-use</i> assays for <ul style="list-style-type: none"> ✓ IL1-β, IFN-γ, IL-2, IL-6, IL-10, IL-4, TNF-α, VEGF, ... ✓ Insulin and glucagon 	<ul style="list-style-type: none"> • Avoids labeling of 1°Abs • Generic pre-labeled 2°Abs (different species available) • Validated for intracellular PTMs, e.g. phosphorylation 	<ul style="list-style-type: none"> • Requires antibody and tracer labeling • Establish competitive antibody binding assays • <i>Ready-to-use</i> assay for hFcRn:Fc protein interactions, e.g. antibodies <ul style="list-style-type: none"> ✓ Lumit™ FcRn Binding Immunoassay



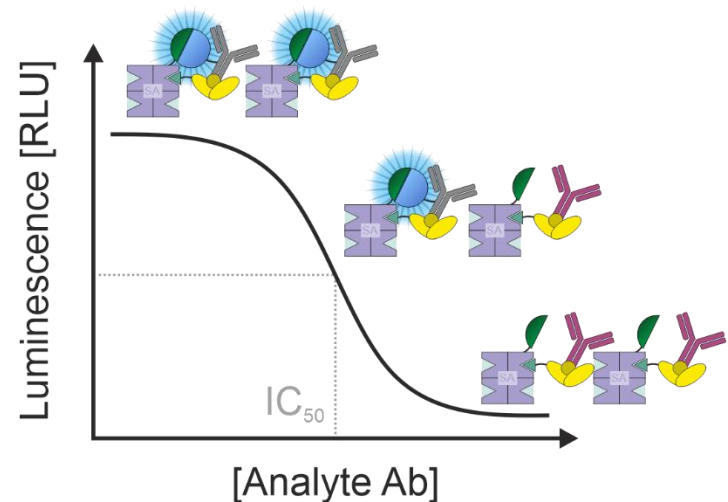
Lumit™ FcRn Binding Immunoassays

Neonatal Fc Receptor Binding Studies without Immobilization Artifacts



FACTS

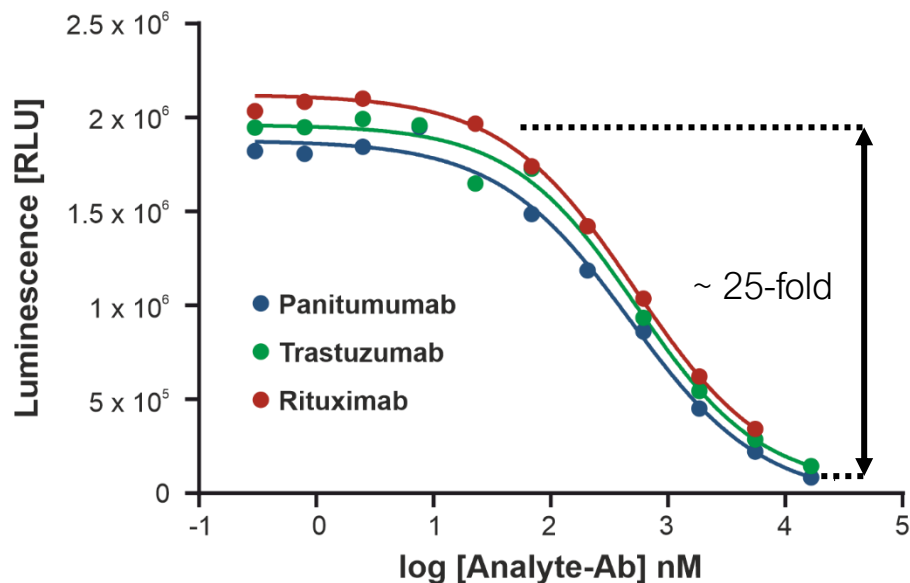
- Tracer: hIgG1 chemically labeled with LgBiT
- Biotinylated hFcRn bound to Streptavidin SmBiT
- Ab-LgBiT:hFcRn SmBiT yields luminescent signal
- Analyte antibodies that bind to hFcRn cause displacement of tracer, i.e. dose dependent decrease in luminescence





Lumit™ FcRn Binding Immunoassays

Analysis of Different Approved Therapeutic Antibodies



Antibody	IC ₅₀ (nM)
Panitumumab	458
Trastuzumab	553
Rituximab	471

OBSERVATION

- Relatively similar IC₅₀ values determined for different therapeutic antibodies as expected
- Good assay window for hIgG:hFcRn binding studies

Summary – Lumit™ Immunoassays

A Bioluminescent Immunodetection Platform based on NanoBiT® Technology

Features of Lumit™ Immunoassays

- Homogenous “add-and read” format can be performed directly on cells
- Fast, 30 to 120 minutes total assay time
- Large dynamic ranges, i.e. no need for sample dilutions
- Use standard plate-reading luminometer
- Scalable to different plate formats (96- and 384-well)
- Amenable to high throughput applications
- Platform includes ready-to-use kits for specific targets as well as reagents to build novel detection assays
- For questions please contact:
erik.bonke@promega.com
- Contact your local sales representative for products & sales relevant information:
www.promega.com/c/local-sales/sales-contacts/

THANK
YOU!