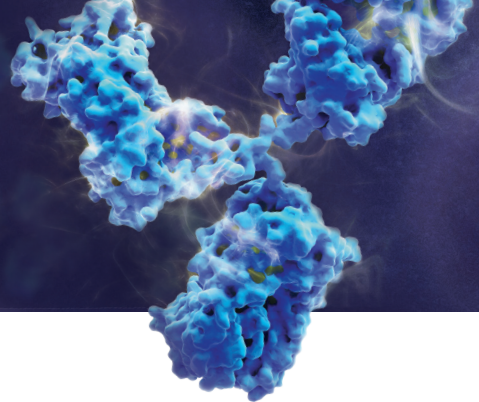


Bioassays for Biologics



Novel Fc Effector and Immune Checkpoint Bioassays for Biologics Drug Discovery and Development

Precise. Accurate. Low variability.

This is a list of Promega bioassay products available through our catalog or through Custom Assay Services (CS numbers).

Fc Effector Bioassays

Bioassay	Cell Components	Catalog or CS#
ADCC Reporter Bioassays		
Complete Kit (WIL2-S) ¹	Effector Cells: FcγRIIIa-V158 Jurkat cells Target Cells: WIL2-S or Raji cells	G7014
Complete Kit (Raji)		G7015
Core Kit ²		G7010
Core Kit (5X)		G7018
Target Kit (WIL2-S) ³		G7013
Target Kit (Raji)		G7016
Propagation Model ⁴		G7102
ADCC Reporter Bioassays, F Variant ⁵		
Core Kit	Effector Cells: FcγRIIIa-F158 Jurkat cells	G9790
Core Kit (5X)		G9798
Propagation Model		G9302
Mouse FcγRIV ADCC Reporter Bioassays		
Complete Kit	Effector Cells: Mouse FcγRIV Jurkat cells	CS1779A05-1
Core Kit		CS1779A01-1
Propagation Model		CS1779A09
FcγRIIIa-H ADCP Reporter Bioassays		
Complete Kit	Effector Cells: FcγRIIIa-H131 Jurkat cells	CS1781A05-1
Core Kit		CS1781A01-1
Propagation Model		CS178101
Anti-TNFα ADCC Bioassays ⁶		
Membrane TNFα Bioassay	Target Cells: Membrane TNFα Cells CHO-K1 cells	CS185502
Membrane TNFα Bioassay, Propagation Model		CS185501
Other Fc Effector Bioassays - Please Enquire		

¹Complete Kits contain Effector Cells, Target Cells, Control Antibody, Cell Culture Medium, Assay Reagents.

²Core Kits contain Effector Cells, Cell Culture Medium, Assay Reagents.

³Target Kits contain Target Cells, Control Antibody.

⁴Propagation Model includes Effector Cells only.

⁵ADCC Reporter Bioassays, F Variant are compatible with the Target Kits listed under "ADCC Reporter Bioassays".

⁶Membrane TNFα CHO-K1 Target Cells are compatible with both the Human and Mouse ADCC Bioassay Effector Cells available in the Core Kits.

Immunotherapy Bioassays

Immune Checkpoint Blockade Bioassays

Bioassay	Cell Components	Catalog or CS#
PD-1/PD-L1 Blockade Bioassays		
PD-1/PD-L1 Blockade Bioassay Kit ¹	PD-1 Effector Cells (Jurkat) PD-L1 aAPC Cells (CHO-K1) ²	CS187111
PD-1/PD-L1 Blockade Bioassay Kit (5X)		CS187119
PD-1/PD-L1 Blockade Bioassay, Propagation Model		CS187109
PD-L1 Negative Cells ³	PD-L1 Negative aAPC Cells (CHO-K1)	CS187104
CTLA-4 Blockade Bioassays		
CTLA-4 Blockade Bioassay	CTLA-4 Effector Cells (Jurkat) aAPC Cells (Raji)	CS186907
CTLA-4 Blockade Bioassay, Propagation Model		CS186906
Other Immune Checkpoint Bioassays - Please Enquire		

¹Kits contain PD-1 Effector Cells, PD-L1 aAPC Cells, Cell Culture Medium, Assay Reagents.

²aAPC (artificial antigen presenting cells) activate the TCR/CD3 complex on Jurkat Effector Cells.

³Negative control cells for the PD-1/PD-L1 Blockade Bioassay.

Immunostimulatory Bioassays

Bioassay	Cell Components	Catalog or CS#
T Cell Activation Bioassays ¹		
T Cell Activation Bioassay (IL-2)	TCR/CD3 Effector Cells (Jurkat)	CS187002
T Cell Activation Bioassay 5X (IL-2)		CS187003
T Cell Activation Bioassay (IL-2), Propagation Model		CS187001
T Cell Activation Bioassay (NFAT)		CS176403
T Cell Activation Bioassay 5X (NFAT)		CS176404
T Cell Activation Bioassay (NFAT), Propagation Model		CS176401
GITR Bioassays		
GITR Bioassay	GITR Effector Cells (Jurkat)	CS184003
CD40 Bioassays		
CD40 Activation Bioassay	CD40 Effector Cells (Ramos)	CS193701
Other Immunostimulatory Receptor Bioassays - Please Enquire		

¹Available with either an IL-2- or NFAT-driven reporter, not both.

Growth Factor and Cytokine Bioassays

Bioassay	Cell Components	Catalog or CS#
TNF α Bioassays		
TNF α Blockade Reporter Bioassay	TNF α Receptor Cells (HEK293)	CS177503
TNF α Blockade Apoptosis Bioassay ¹	TNF α Receptor Cells (U937)	CS1324A05
Membrane TNF α Bioassay	Target Cells: Membrane TNF α Cells CHO-K1 cells	CS185502
Membrane TNF α Bioassay, Propagation Model		CS185501
VEGF Bioassays		
VEGF Blockade Bioassay	KDR Cells (HEK293)	CS181402
Other Growth Factor and Cytokine Receptor Bioassays - Please Enquire		

¹Designed for use with Promega's Caspase-Glo® 3/7 Assay Reagents (Cat.# [G8091](#)).

For more information on bioassays, visit: www.promega.com/BioAssay



Build It, Buy It or Customize Your Assay

Our Custom Assay Services Team can build a complete biology-driven, technology-enabled custom solution.

For more information on custom assays visit www.promega.com/CAS or email CAS@promega.com