

Products + Services



Quantus™ Fluorometer

A dual-channel fluorometer for quantifying nucleic acids. This compact instrument is simple to operate, has preprogrammed settings for Promega QuantiFluor® Dye Systems and allows users the flexibility to create their own methods and quantitation settings for other dyes.

[Learn more](#)

[Product Details + Ordering](#)

ALL CATEGORIES

Biobanking

DNA and RNA Quantitation, DNA Extraction for Biobanks, Sample ID and Mixed Sample Detection

Biochemicals and Labware

Biochemical Buffers and Reagents, Nucleic Acids, Nucleic Acids Purification Accessories, Tips and Accessories

Biologics

Bioanalytical Tools, Bioassays for Biologics, Protein Manipulation Tools

Cell Health and Metabolism

ADME Assays, Apoptosis Assays, Cell Viability Assays, Cytotoxicity Assays, View All

Cell Signaling

AMP Detection System, GPCR Assays, Growth Factors, Histone Deacetylase Assays, View All

Cloning and DNA Markers

Cloning Tools and Competent Cells, Molecular Weight Markers, PCR Cloning, Restriction Enzymes

DNA and RNA Purification

Genomic DNA Purification Kits, Plasmid Purification, RNA Purification, Nucleic Acids Purification Accessories, View All

Drug Discovery

ADME Assays, Apoptosis Assays, Cell Viability Assays, Custom Assay Materials, View All

Epigenetics

Cell-Based and Biochemical Assays, DNA Methylation Analysis, Protein Analysis and Complex Purification

Genetic Identity

Automation for Genetic Identity, DNA Quantitation for Genetic Identity, Sample Preparation for Genetic Identity, STR Analysis for Forensic and Paternity Testing

Imaging and Immunological Detection

Cellular Imaging with HaloTag®, ELISAs and Antibodies, In vivo Imaging

Industrial and Environmental Monitoring

DNA Purification From Food, Microbial Detection and Quantitation, Protein Deamidation Detection

Instruments

Fluorometers, HSM 2.0 Instrument, Luminometers, Maxwell® Research Systems, View All

Molecular Diagnostics

Maxwell® Research Systems, Maxwell® CSC System for IVD Use, Systems for Molecular Diagnostics Research, Reagents for Molecular Diagnostics Labs

PCR

Hot-Start PCR, Routine PCR, Long PCR, qPCR and RT-qPCR, View All

Protein Expression and Mass Spectrometry

Eukaryotic Cell-Free Protein Expression, Prokaryotic Cell-Free Protein Expression, Cell-Based Protein Expression, Glycosidases, View All

Protein Purification and Interactions

HaloTag® Protein Purification, GST Protein Purification, Magnetic Systems for Purification of Antibodies and Affinity-Tagged Proteins, Protein Interactions, View All

Reporter Assays and Transfection

In vivo Imaging, Reporter Assays, Reporter Vectors and Cell Lines, Transfection Reagents

RNA Analysis

In Vitro Transcription, RNA Interference, RNA Purification, RT-PCR

Stem Cell Research

Cell Line Authentication, Stem Cell Functional Analysis, Stem Cell Health, Stem Cell Sample Preparation, View All

STR Analysis

Sample ID and Mixed Sample Detection, Cell Line Authentication, STR Analysis for Forensic and Paternity Testing

Vectors

Bacterial Expression Vectors, Cell-Free Expression Vectors, Mammalian Expression Vectors, Reporter Vectors, View All