

# NEW MSD-ECL Protein Detection Service

**Validated Immunoassays for Reproducible Results. Unbeatable Matrix Tolerance for a Range of Applications.**

The recent addition of the MESO SCALE DISCOVERY® Electrochemiluminescence (MSD-ECL) platform to PBL Assay Science's existing service offerings enhances our capability to provide broad analyte quantification in normal and disease samples in singleplex and multiplex formats, and greatly expands the menu of crucial analytes.

## Broad Range of Applications:

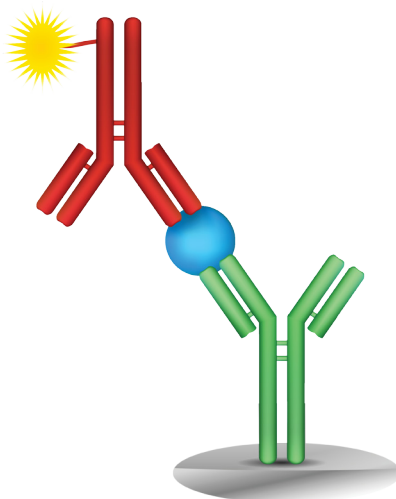
- Cytokines & Chemokines
- Cardiac Biomarkers
- Immunogenicity
- Immunology
- Inflammation
- Intracellular Signaling
- Metabolic
- Neurodegeneration
- Oncology & Cancer
- Toxicology
- ...and more

Multiplex immunoassays on the MSD® platform maintain the sensitivity and performance offered by singleplex ELISAs while providing additional benefits such as cost-savings and the potential to target several analytes in a single sample. The ability to quantify up to 10 compatible analytes while requiring no more than 25 µl of neat sample ensures efficient use of precious sample. Entrusting your sample testing analysis to PBL's experienced team of dedicated assay services scientists and quality control experts ensures trustworthy and accurate results.

PBL offers high precision protein quantification services for the detection of biomarkers in a broad range of sample matrices, including, but not limited to, serum, plasma, and tissue culture media for Human, Non-Human Primate (NHP), Mouse, Rat, and Canine samples. Inquire with our Assay Services team to learn more about our capabilities and how we can help advance your research with the MSD platform.

## MSD-Electrochemiluminescence (ECL) Technology Overview

Unique detection technology provided by the MSD platform utilizes SULFO-TAG™ labels which emit light upon electrochemical stimulation. The intensity of light generated is captured via CCD camera and pixel intensities quantified. This system provides the basis for achieving measurement of a several log range of biomarker expression levels in a variety of matrices.



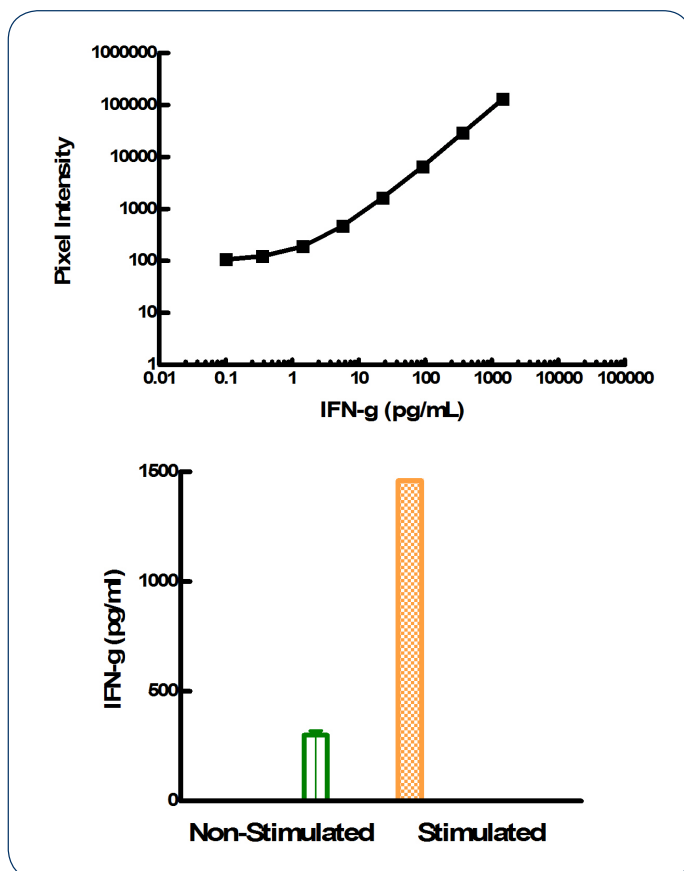
# MSD® Electrochemiluminescence (ECL)

High Performance Protein Detection Services in Singleplex and Multiplex Formats

**Broad Dynamic Range. Minimal Background. High Sensitivity.**

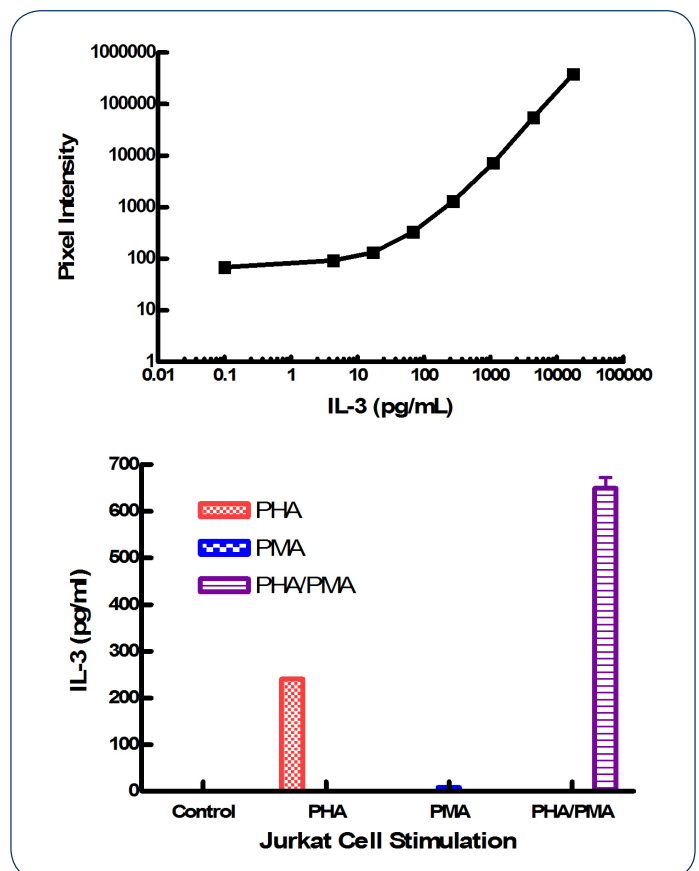
**Enhanced signal capture achieved through luminescent detection.**

- Ability to target low analyte levels in **small sample volumes** empowered by ECL technology
- Robust **matrix tolerance** including serum, plasma, and TCM, as well as various complex biological matrices
- Obtain accurate measurement of normal and elevated analyte within the same sample with **5-Log+ dynamic range**
- Patterned arrays and parallel biological processing offer optimal **multiplex** functionality
- Over 180 **validated** fit-for-purpose singleplex and preconfigured multiplex panel ELISAs available with guaranteed performance specifications and lot-to-lot consistency for reproducible results



(top) Human IFN-g standard curve exhibits a wide dynamic range allowing for sensitive measurement in undiluted samples.

(bottom) Human IFN-g released from NK-92 cells upon exposure to IL-2.



(top) Human IL-3 standard curve demonstrates the broad dynamic range characteristic of MSD immunoassays.

(bottom) IL-3 production by Jurkat cell line after stimulation by PHA, PMA, and both agents.