

InCELL Hunter™ A549 CSF1R Kinase Cell Line

Catalog Number: 96-0047C15 **Lot Number:** See Vial
Contents: 2 vials, 1 x 10⁶ cells per vial in 1 mL

Background

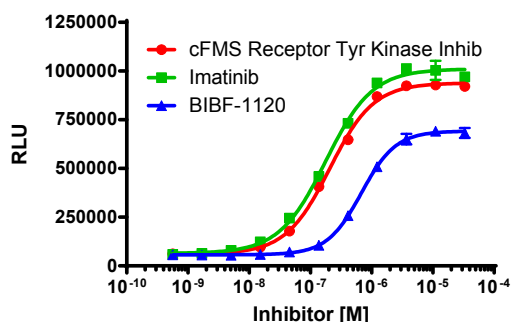
The InCELL Hunter technology measures intracellular compound-target engagement by detecting changes in protein stability due to compound binding. In this assay, the target is fused to the enhanced ProLabel® (ePL) enzyme fragment and expressed in the selected cell background. The amount of fusion is quantified through the addition of substrate and complementary enzyme acceptor fragment (EA). The small size, but robust activity of the ePL tag makes it an ideal fusion partner. The small disordered peptide has minimal effects on target half-life, and generates high levels of enzyme activity, resulting in a sensitive measurement of protein levels.

Product Information

Target: CSF1R Amino Acid Range: (aa 560 - 937)
Description: colony-stimulating factor 1 receptor
Accession Number: NM_005211.3
Target Species: Human
ProLabel™ Tag: ePL **ProLabel™ Tag located at (terminus):** C-Terminus
Cell Type: A549
Storage: Short term (<24 h): Store at -80°C; Long term (>24 h): Store in vapor phase of liquid nitrogen.

Functional Performance

Cells were seeded in a 384-well plate, incubated at 37°C and 5% CO₂ followed by stimulation with a control compound, as defined in the assay conditions below. Following treatment, signal was detected using the InCELL Hunter Detection Kit according to the recommended protocol using the conditions specified below. Please refer to page 2 for recommended assay reagents, detection reagents, and control compounds.



Cell Number/Well: 5000
Cell Seeding Time (Hours): 24
Control Compound: cFMS Receptor Tyrosine Kinase Inhibitor
Compound Incubation Time (minutes): 360
Compound Incubation Temperature (°C): 37
EC₅₀ of Compound (nM): 207.0
Signal:Background at Compound E_{max}: 15.5
1:4 EA Dilution Required? Yes
Detection Incubation Time (hours): 1

Passage Stability

This cell line has been confirmed to be stable through 10 passages with no significant drop in assay window or change in EC₅₀.

Mycoplasma Testing

This lot was tested and found to be free of mycoplasma contamination. Data available upon request.

Required Materials

The following additional materials are required but not provided:

Product Use*	Product Description	Catalog Number
Detection	InCELL Detection Kit	96-0079
Cell Culture	AssayComplete™ Cell Culture Kit-105	92-3105G
Cell Plating	AssayComplete™ Cell Plating 0 Reagent	93-0563R0A
Cell Detachment	AssayComplete™ Cell Detachment Reagent	92-0009
Cell Thawing	AssayComplete™ Thawing Reagent T1	92-4101TR
Cell Freezing	AssayComplete™ Freezing Reagent F1	92-5101FR

*Please inquire about our cell line-specific AssayComplete Starter Packs to get you started with your cell culture needs.

Required Antibiotics

Antibiotic Name	Concentration (µg/mL)	Catalog Number
AssayComplete™ Puromycin	Not Applicable	Not Applicable
AssayComplete™ Hygromycin B	Not Applicable	Not Applicable
AssayComplete™ G418	500	92-0030

Additional Compound Information

Control Inhibitor: cFMS Receptor Tyrosine Kinase Inhibitor

Vendor: DiscoverRx® (Catalog No. 92-1240)

For order placement or technical support, please call 1.866.448.4864 (North America) or +44.121.260.6142 (Europe) or e-mail info@discoverx.com. For additional information, please visit discoverx.com.

Limited Use License Agreement

These products may be covered by issued US and/or foreign patents, patent application and subject to Limited Use Label License.

Please visit discoverx.com/license for a list of products that are governed by limited use label license terms and relevant patent and trademark information.