

PathHunter® Jurkat ICOS Signaling Assay

Materials Provided

Catalog Number: 93-1127C19

Lot Number: See Vial

Background

Checkpoint Signaling assays are engineered to co-express a ProLink™ (PK) tagged immune checkpoint receptor and an Enzyme Acceptor (EA) tagged SH2 domain. Ligand engagement leads to receptor activation and phosphorylation, resulting in SH2-EA recruitment to the receptor, and forcing complementation of the two β-galactosidase enzyme fragments (EA and PK). The resulting functional enzyme hydrolyzes substrate to generate a chemiluminescent signal.

Product Information

Target Receptor: ICOS

Target Accession #: NM_012092.3

Description: Inducible T-cell costimulator

Species: Human

SH2 Domain: P85

SH2 Accession #: NM_181523.2

Amino Acid Range: 320 - 435

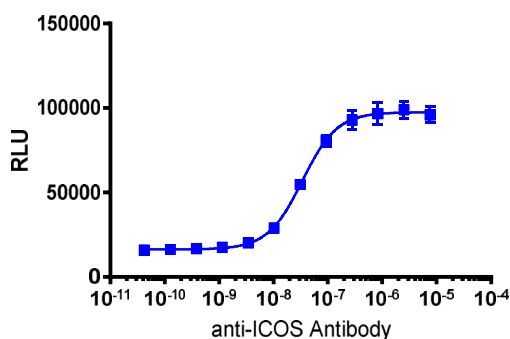
Cell Type: Jurkat (Suspension)

Contents: 4 x 10⁶ cells per vial in 1 mL

Storage: Short term (<24 h): Store at -80°C; Long term (>24 h): Store in vapor phase of liquid nitrogen.

Functional Performance

Signaling assay cells were seeded in a 96 well plate and incubated at 37°C and 5% CO₂ for the indicated time specified below. Following agonist incubation, signal was detected using the PathHunter Detection Kit according to the recommended protocol. Please refer to page 2 for recommended assay reagents, detection reagents, and control compounds.?



Signaling Cell Number/Well:	30000
Signaling Cell Seeding Time (hours):	4
Control Agonist:	Anti-ICOS antibody
Agonist Incubation Time (minutes):	120
Agonist Incubation Temperature (°C):	RT
EC₅₀ for compound stimulation (ng/mL):	34.3
Signal:Background at agonist E_{max}:	6.3

Detection Incubation Time (hrs.)	1
---	---

Passage Stability

This cell line has been confirmed to be stable through a minimum of 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	PathHunter® Bioassay Detection Kit	93-0933
Cell Culture	AssayComplete™ Cell Culture Kit-101	92-3101G
Cell Plating	AssayComplete™ Cell Plating 0 Reagent	93-0563R0A
Cell Detachment	Not Applicable	Not Applicable
Cell Thawing	AssayComplete™ Thawing Reagent T6	92-4106TR
Cell Freezing	AssayComplete™ Freezing Reagent F5	92-5105FR

*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	250	92-0029
AssayComplete™ G418	500	92-0030

Additional Ligand Information

Control Agonist: Anti-ICOS antibody

Vendor: Eurofins DiscoverX® (Catalog No. 92-1339)

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 <https://www.discoverx.com/legal-policies/> for limited-use label license agreement and trademark information for this product.