

Human c-MET/HGFR Reporter Assay System

This is an all-inclusive cell-based luciferase reporter assay kit targeting the the Human c-MET / Hepatocyte Growth Factor Receptor (c-MET/HGFR). INDIGO's c-MET/HGFR reporter assay utilizes proprietary mammalian cells that have been engineered to provide constitutive expression of the OXTR. In addition to c-MET/HGFR Reporter Cells, this kit provides two optimized media for use during cell culture and in diluting the user's test samples, a reference agonist, Luciferase Detection Reagent, and a cell culture-ready assay plate. The principal application of this assay is in the screening of test samples to quantify any functional activity, either agonist or antagonist, that they may exert against c-MET/HGFR.

Ordering Information

[Order Online »](#)

NGO-IB30301-0096	96 wells
NGO-IB30302-0384	384 wells

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name c-MET / Hepatocyte Growth Factor Receptor

Contents

- Human c-MET/HGFR Reporter Cells
- Cell Recovery Medium (CRM)
- Compound Screening Medium (CSM)
- HGF, (ref. agonist; in PBS/0.1%BSA)
- Detection Substrate
- Detection Buffer
- White, sterile, cell-culture ready assay plate

Species Reactivity Human



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com