

# Human G-CSFR Reporter Assay System

This is an all-inclusive cell-based luciferase reporter assay kit targeting the the Human Granulocyte Colony-Stimulating Factor Receptor (G-CSFR). INDIGO's G-CSFR reporter assay utilizes proprietary mammalian cells that have been engineered to provide constitutive expression of the G-CSFR. In addition to G-CSFR 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 G-CSFR.

## Ordering Information

[Order Online »](#)

NGO-IB30101-0096	96 wells
NGO-IB30102-0384	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** Granulocyte Colony-Stimulating Factor Receptor

**Contents**

- Human G-CSFR Reporter Cells
- Cell Recovery Medium (CRM)
- Compound Screening Medium (CSM)
- G-CSF, (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](mailto:info-usa@enzolifesciences.com)

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

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

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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