

Human GM-CSR Reporter Assay System

Ordering Information

[Order Online »](#)

NGO-IB30001-0096	96 wells
NGO-IB30002-0384	384 wells

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Granulocyte-Macrophage Colony-Stimulating Factor Receptor

Contents This Granulocyte-Macrophage Colony-Stimulating Factor Receptor (GM-CSFR) assay kit is an all-inclusive luciferase reporter assay system that includes, in addition to GM-CSFR Reporter Cells, optimized culture medium for use in reviving the cryopreserved cells, a culture medium for use in preparing test sample treatments, the physiological activator GM-CSF as a positive control, Luciferase Detection Reagents, a cell culture-ready assay plate, and a detailed protocol.



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