

Human EGFR1 Reporter Assay System

Ordering Information

[Order Online »](#)

NGO-IB13001-0096	96 wells
NGO-IB13001-0032	3x32 assays
NGO-IB13002-0384	384 wells

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Shipping

Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name

Epidermal Growth Factor Receptor 1

Contents

INDIGO's EGFR1 assay kit provides the convenience of an all-inclusive cell-based assay system. In addition the the EGFR1 Reporter Cells, provided are two optimized media for use in recovering the cryopreserved cells and for diluting test samples. Also included is the reference agonist EGF, Luciferase Detection Reagents, and a cell culture-ready assay plate.



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