

Rat GR Reporter Assay System

INDIGO's Rat Glucocorticoid Receptor (nr3c1) nuclear receptor assay utilizes proprietary non-human cells engineered to provide constitutive high-level expression of full-length, unmodified Rat Glucocorticoid Receptor (nr3c1), a ligand-dependent transcription factor, commonly referred to as GR.

Ordering Information

[Order Online »](#)

NGO-R00201-0096	96 wells
NGO-R00201-0032	3x32 assays

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Glucocorticoid Receptor

Contents This kit product is an all-inclusive assay system that includes, in addition to rGR Reporter Cells, two optimized media for use during cell culture and (optionally) in diluting the test samples, a reference agonist, Luciferase Detection Reagent, 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