

Zebrafish GR Reporter Assay System

INDIGO's Zebrafish Glucocorticoid Receptor (nr3c1) nuclear receptor assay utilizes proprietary mammalian cells engineered to provide constitutive highlevel expression of full-length, unmodified Zebrafish (Danio rerio) Glucocorticoid Receptor (nr3c1), a ligand-dependent transcription factor, commonly referred to as GR.

Ordering Information

[Order Online »](#)

NGO-Z00201-0096	96 wells
NGO-Z00201-0032	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

INDIGO Bioscience's receptor assays are all-inclusive cell-based assay systems. In addition to zGR Reporter Cells, this kit provides two optimized media for use during cell culture and in diluting the user's test samples, the reference agonist Dexamethasone, Luciferase Detection Reagent, 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