

Zebrafish TR β Reporter Assay System

INDIGO's Zebrafish Thyroid Hormone Receptor Beta Reporter Assay System utilizes proprietary human cells engineered to provide constitutive, high-level expression of Zebrafish (*Danio rerio*) Thyroid Hormone Receptor Beta (nr1a2), a ligand-dependent transcription factor.

INDIGO's Reporter Cells include the luciferase reporter gene functionally linked to a TR β -responsive promoter. Thus, quantifying changes in luciferase expression in the test sample-treated reporter cells provides a sensitive surrogate measure of changes in zTR β activity. The principal application of this assay is in the screening of test samples to quantify any functional bioactivity that they may exert against zebrafish AR. In particular, zebrafish reporter assays are frequently used in the monitoring of environmental samples for the presence of biohazardous chemical pollutants, such as endocrine disruptors.

Ordering Information

[Order Online »](#)

NGO-Z01101-0032	3x32 assays
NGO-Z01101-0096	96 wells

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Thyroid Hormone Receptor β

Contents This kit product is an all-inclusive assay system that includes, in addition to zTR β 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