

Zebrafish PPAR γ Reporter Assay System

INDIGO's Zebrafish Peroxisome Proliferator Activated Receptor Gamma Reporter Assay System utilizes proprietary mammalian cells engineered to provide constitutive, high-level expression of Zebrafish (*Danio rerio*) PPAR γ (nr1c3), a ligand-dependent transcription factor.

Ordering Information

[Order Online »](#)

NGO-Z00101-0096	96 wells
NGO-Z00101-0032	3x32 assays

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Shipping

Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name

Peroxisome Proliferator-Activated Receptor Gamma

Contents

This kit product is an all-inclusive assay system that includes, in addition to zPPAR γ 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