

Cyn Monkey PPAR α Reporter Assay System

INDIGO's Cynomolgus Monkey Peroxisome Proliferator-Activated Receptor Alpha (cPPAR α) Reporter Assay System utilizes proprietary mammalian cells engineered to provide constitutive, high-level expression of Cyn Monkey PPAR α (nr1c1), a ligand-dependent transcription factor. Because these cells incorporate a responsive luciferase reporter gene, quantifying expressed luciferase activity provides a sensitive surrogate measure of cPPAR α activity in treated cells. The primary application of this reporter assay system is in the screening of test samples to quantify functional activity, either agonist or antagonist, that they may exert against cynomolgus monkey PPAR α .

Ordering Information

[Order Online »](#)

NGO-C00111-0032	3x32 assays
NGO-C00111-0096	96 wells

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Shipping

Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name

Peroxisome Proliferator-Axivated Receptor α

Contents

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