

# Cyn Monkey PPAR $\gamma$ Reporter Assay System

INDIGO's Cynomolgous Monkey PPAR gamma (cPPAR $\gamma$ ; nr1c3) Assay utilizes proprietary non-human mammalian cells engineered to provide high-level expression of cPPAR $\gamma$ . Reporter Cells also incorporate an cPPAR $\gamma$ -responsive luciferase reporter gene, therefore, quantifying expressed luciferase activity provides a sensitive surrogate measure of cPPAR $\gamma$  activity in the treated cells. The principle application of this reporter assay system is in the screening of test samples to quantify any functional activity, either agonist or antagonist, that they may exert against cPPAR $\gamma$ .

## Ordering Information

[Order Online »](#)

NGO-C00101-0032	3x32 assays
NGO-C00101-0096	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  $\gamma$

**Contents**

This kit product is an all-inclusive assay system that includes, in addition to cPPAR $\alpha$  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](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)