

ROR γ (h) Reporter Assay System

Human ROR γ compound screening assay

This nuclear receptor assay system utilizes proprietary non-human mammalian cells engineered to provide high-level expression of a hybrid form of the Human RAR-related Orphan Receptor Gamma. The N-terminal DNA binding domains (DBD) of the native ROR γ and ROR γ t receptors have been substituted with that of the yeast GAL4-DBD. Hence, the GAL4-ROR γ hybrid receptor expressed in these reporter cells will not discern any functional differences that may exist between the native isoform 1 and isoform 2 receptors. The principle application of this reporter assay system is in the screening of test samples to quantify inverse-agonist or antagonist activities that they may exert against human ROR γ .

Ordering Information

[Order Online »](#)

NGO-IB04001-0096	96 wells
NGO-IB04001-0032	3x32 assays
NGO-IB04002-0384	384 wells

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Human RAR-related Orphan Receptor Gamma Assay System, NR1F3, RORC, RORg

Application HTS

Application Notes The principle application of this reporter assay system is in the screening of test samples to quantify inverse-agonist or antagonist activities that they may exert against human ROR γ .

Contents

- ROR γ Reporter Cells
- Cell Recovery Medium (CRM)
- Compound Screening Medium (CSM)
- Ursolic Acid (reference inverse-agonist for RORg), 10 mM (in DMSO)
- Detection Substrate
- Detection Buffer
- 96-well assay plate (white, sterile, cell-culture ready) (96 assays)
- 96-well plate frame + Snap-in 8-well strips (3 x 32 assays)
- 384-well assay plate (384 assays)



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