

# RAR $\beta$ (h) Reporter Assay System

## Human RAR $\beta$ compound screening assay

This nuclear receptor assay system utilizes proprietary non-human mammalian cells engineered to provide constitutive, high-level expression of Human Retinoic Acid Receptor, Beta (NR1B2), a ligand-dependent transcription factor commonly referred to as RARB or RAR $\beta$ . Because these cells incorporate a responsive luciferase reporter gene, quantifying expressed luciferase activity provides a sensitive surrogate measure of RAR $\beta$  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 human RAR $\beta$ .

### Ordering Information

[Order Online »](#)

NGO-IB02101-0096	96 wells
NGO-IB02101-0032	3x32 assays
NGO-IB02102-0384	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 Retinoic Acid Receptor, Beta (RAR $\beta$ ) Reporter Assay System, NR1B2, RARB, RAR $\beta$

**Application** HTS

**Application Notes** 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 human RAR $\beta$ .

**Contents**

- RAR $\beta$  Reporter Cells
- Cell Recovery Medium (CRM)
- Compound Screening Medium (CSM)
- trans-Retinoic Acid (reference agonist for RAR $\beta$ ), 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](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)