

# Human ADRB1 Reporter Assay System

This is an all-inclusive cell-based luciferase reporter assay kit targeting the Human Adrenoreceptor Beta 1 (ADRB1). INDIGO's ADRB1 reporter assay utilizes proprietary mammalian cells that have been engineered to provide constitutive expression of the ADRB1. In addition to ADRB1 Reporter Cells, this kit provides two optimized media for use during cell culture and in diluting the user's test samples, a reference agonist, Luciferase Detection Reagent, and a cell culture-ready assay plate. The principal application of this assay is in the screening of test samples to quantify any functional activity, either agonist or antagonist, that they may exert against ADRB1.

## Ordering Information

[Order Online »](#)

NGO-IB32001-0096	96 wells
NGO-IB32002-0384	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              Adrenoreceptor Beta 1

Contents

- Human ADRB1 Reporter Cells
- Cell Recovery Medium (CRM)
- Compound Screening Medium (CSM)
- Norepinephrine, (ref. agonist; in DMSO)
- Detection Substrate
- Detection Buffer
- White, sterile, cell-culture ready assay plate

Species Reactivity              Human



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)