

Human ADRA2A Reporter Assay System

This is an all-inclusive cell-based luciferase reporter assay kit targeting the Human Adrenoceptor Alpha 2A (ADRA2A). INDIGO's ADRA2A reporter assay utilizes proprietary mammalian cells that have been engineered to provide constitutive expression of the ADRA2A. In addition to ADRA2A 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 ADRA2A.

Ordering Information

[Order Online »](#)

NGO-IB34001-0096	96 wells
NGO-IB34002-0384	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 Adrenoceptor Alpha 2A

Contents

- Human ADRA2A Reporter Cells
- Cell Recovery Medium (CRM)
- Compound Screening Medium (CSM)
- Brimonidine, (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

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