

Human AVPR1B Reporter Assay System

This is an all-inclusive cell-based luciferase reporter assay kit targeting the Human Arginine Vasopressin Receptor 1B (AVPR1B). INDIGO's AVPR1B reporter assay utilizes proprietary mammalian cells that have been engineered to provide constitutive expression of the AVPR1B.

Ordering Information

[Order Online »](#)

NGO-IB37101-0096	96 wells
NGO-IB37102-0384	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 Arginine Vasopressin Receptor 1B

Contents

- Human AVPR1B Reporter Cells
- Cell Recovery Medium (CRM)
- Compound Screening Medium (CSM)
- Vasopressin, (ref. agonist; in PBS+0.1%BSA)
- Detection Substrate
- Detection Buffer
- White, sterile, cell-culture ready assay plate

Species Reactivity Human



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info@enzolifesciences.com
usa@enzolifesciences.com

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info@enzolifesciences.com
be@enzolifesciences.com

France
Phone: +33 472 440 655
info@enzolifesciences.com
fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info@enzolifesciences.com
de@enzolifesciences.com

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