

# Human AVPR2 Reporter Assay System

This is an all-inclusive cell-based luciferase reporter assay kit targeting the Human Arginine Vasopressin Receptor 2 (AVPR2). INDIGO's AVPR1B reporter assay utilizes proprietary mammalian cells that have been engineered to provide constitutive expression of the AVPR2.

## Ordering Information

[Order Online »](#)

NGO-IB37201-0096	96 wells
NGO-IB37202-0384	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 2

**Contents**

- Human AVPR2 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](mailto:info@enzolifesciences.com)  
[usa@enzolifesciences.com](mailto:usa@enzolifesciences.com)

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

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info@enzolifesciences.com](mailto:info@enzolifesciences.com)  
[be@enzolifesciences.com](mailto:be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info@enzolifesciences.com](mailto:info@enzolifesciences.com)  
[fr@enzolifesciences.com](mailto:fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info@enzolifesciences.com](mailto:info@enzolifesciences.com)  
[de@enzolifesciences.com](mailto:de@enzolifesciences.com)

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