

Human CB1R Reporter Assay System

This assay uses proprietary human cells that provide constitutive expression of the Human Cannabinoid Type 1 Receptor, also referred to as CB1R.

Ordering Information

[Order Online »](#)

NGO-IB19001-0032	3x32 assays
NGO-IB19001-0096	96 wells
NGO-IB19002-0384	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 Cannabinoid Type 1 Receptor

Contents This Human Cannabinoid Type 1 Receptor (CB1R) assay kit is an all-inclusive assay system that includes addition to CB1R Reporter Cells, two optimized media for use in recovering the cryopreserved cells and for diluting test samples, the reference activator synthetic cannabinoid CP-55,940, Luciferase Detection Reagents, and a cell culture-ready assay plate.



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