

# Human VEGFR2 Reporter Assay System

This Human type II Vascular Endothelial Growth Factor Receptor (VEGFR2) kit product is an all-inclusive assay system that includes addition to VEGFR2 Reporter Cells, two optimized media for use in recovering the cryopreserved cells and for diluting test samples, the reference activator human VEGF-A (165 aa isoform, non-glycosylated), Luciferase Detection Reagents, and a cell culture-ready assay plate.

## Ordering Information

[Order Online »](#)

NGO-IB15001-0096	96 wells
NGO-IB15001-0032	3x32 assays
NGO-IB15002-0384	384 wells

## Manuals, SDS & CofA

[View Online »](#)

# Handling & Storage

Shipping

Dry Ice

**Regulatory Status** RUO - Research Use Only

# Product Details

**Alternative Name**

Vascular endothelial growth factor receptor 2



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)