

# Human IGF-1R Reporter Assay System

## Ordering Information

[Order Online »](#)

NGO-IB29001-0096	96 wells
NGO-IB29002-0384	384 wells

## Manuals, SDS & CofA

[View Online »](#)

# Handling & Storage

Shipping

Dry Ice

# Regulatory Status

RUO - Research Use Only

# Product Details

Alternative Name

Insulin-Like Growth Factor Receptor 1

Contents

This Insulin-Like Growth Factor Receptor 1 (IGF-1R) assay kit is an all-inclusive luciferase reporter assay system that includes, in addition to IGF-1R Reporter Cells, optimized culture media for use in recovering the cryopreserved cells a culture medium for use in preparing test sample treatments, the physiological activator Insulin as a positive control, Luciferase Detection Reagents, a cell culture-ready assay plate, and a detailed protocol.



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)