

# Human FGFR1/2 Reporter Assay System

INDIGO's assay utilizes proprietary human cells that have been engineered to provide constitutive expression of the Human Fibroblast Growth Factors 1 and 2 (FGFR1/2). FGFR1 and FGFR2 are both single-pass transmembrane receptors that contain respective extracellular ligand-binding domains, transmembrane domains, and intracellular tyrosine kinase domains.

## Ordering Information

[Order Online »](#)

NGO-IB21001-0032	3x32 assays
NGO-IB21001-0096	96 wells
NGO-IB21002-0384	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**              Fibroblast Growth Factor Receptors 1 & 2

**Contents**                      This Human Fibroblast Growth Factor Receptor 1 and 2 (FGFR1/2) assay kit is an all-inclusive assay system that includes addition to FGFR1/2 Reporter Cells, an optimized culture medium for use in reviving the cryopreserved cells, a culture medium supplemented with heparin for use in preparing test sample treatments, the physiological activator FGF-Acidic (aka FGF1), Luciferase Detection Reagents, and a cell culture-ready assay plate.



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)