

# Human FGFR4 Reporter Assay System

This is an all-inclusive cell-based luciferase reporter assay kit targeting the Human Fibroblast Growth Factor Receptor 4 (FGFR4). INDIGO's FGFR4 reporter assay utilizes proprietary mammalian cells that have been engineered to provide constitutive expression of the Human Fibroblast Growth Factor Receptor 4. In addition to FGFR4 Reporter Cells, this kit provides two optimized media for use during cell culture and in diluting the user's test samples, a reference agonist, Luciferase Detection Reagent, and a cell culture-ready assay plate. The principal application of this assay is in the screening of test samples to quantify any functional activity, either agonist or antagonist, that they may exert against human FGFR4. This kit provides researchers with clear, reproducible results, exceptional cell viability post-thaw, and consistent results lot to lot.

The family of Fibroblast Growth Factors (FGFs) comprise approximately 23 members that are related by core sequence and structure conservation, with the majority of FGFs being secreted signaling proteins. Secreted FGFs are predominantly autocrine and paracrine factors, with only three members evolved to function as endocrine factors.

FGFs bind and activate four FGF Receptors (FGFR1-4) which, themselves, are members of the family of high-affinity tyrosine kinase receptors. Heparin and heparin sulfate proteoglycans (HSPGs) are essential cofactors for paracrine FGF (e.g., FGF-1 and FGF-2) interactions with FGFRs. The association between paracrine FGFs and HSPGs ensures their limited diffusion and enhanced FGFR binding specificity.

## Ordering Information

[Order Online »](#)

NGO-IB43001-0096	96 wells
NGO-IB43002-0384	384 wells

## Manuals, SDS & CofA

[View Online »](#)

- Ready to Use Upon Receipt
- Includes All Needed Components
- Contains Transfected Reporter Cells
- Eliminates Cell Licensing Fees
- Clear, Reproducible Results
- Consistent Results Lot to Lot

## Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Fibroblast Growth Factor Receptor 4

**Contents**

- FGFR4 Reporter Cells
- Cell Recovery Medium (CRM)
- Compound Screening Medium (CSM)
- FGF-2, (ref. activator; 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)