

FGFR/ β -Klotho Reporter Assay System

INDIGO's Human Fibroblast Growth Factor Receptor Receptor 1c and β -Klotho (FGFR/ β -Klotho) assay utilizes proprietary human cells that have been engineered to provide constitutive expression of the Human Fibroblast Growth Factor Receptor 1c and β -Klotho. FGFR1c and β -Klotho are both single-pass transmembrane proteins. FGFR1c has an extracellular ligand-binding domain, transmembrane domain, and intracellular tyrosine kinase domain. It has been established that FGFR association with the co-receptor β -Klotho generates a scaffold that is essential for endocrine growth factor binding interactions, such as those with FGF-21 and FGF-19. Following growth factor binding, the activated tyrosine kinase activities of the FGFR initiate intracellular signaling cascades that may include RAS-MAPK, PI3-AKT, PLC γ and/or STAT pathways. It is FGFR1c/ β -Klotho signal transduction via the RAS-MAPK-ERK1/2 cascade that is exploited by the reporter cells in this assay.

Ordering Information

[Order Online »](#)

NGO-IB22001-0032	3x32 assays
NGO-IB22001-0096	96 wells
NGO-IB22002-0384	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 Receptor 1c and β -Klotho

Application Notes The principal application of this reporter assay is in the screening of test materials to quantify any functional interactions, either activating or inhibitory, that they may exert against FGFR1c/ β -Klotho or the coupled RAS-MAPK pathway.

Contents This Human Fibroblast Growth Factor Receptor 1c and β -Klotho (FGFR/ β -Klotho) assay kit is an all-inclusive assay system that includes addition to FGFR/ β -Klotho Reporter Cells, an optimized culture medium for use in reviving the cryopreserved cells, a culture medium for use in preparing test sample treatments, the physiological endocrine activator FGF-21, 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