

Human FGFR1/ α -Klotho Reporter Assay System

This is an all-inclusive cell-based luciferase reporter assay kit targeting the Human Fibroblast Growth Factor Receptor 1c and α -Klotho (FGFR/ α -Klotho). INDIGO's FGFR/ α -Klotho reporter assay utilizes proprietary mammalian cells that have been engineered to provide constitutive expression of the Human Fibroblast Growth Factor Receptor 1c and α -Klotho.

Ordering Information

[Order Online »](#)

NGO-IB40001-0096	96 wells
NGO-IB40002-0384	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 1/ α -Klotho

Contents

- FGFR/ α -Klotho Reporter Cells
- Cell Recovery Medium (CRM)
- Compound Screening Medium (CSM)
- FGF-23, (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-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