

Human ERRb Reporter Assay System

This Estrogen-related Receptor Beta (ERR β) assay kit is an all-inclusive firefly luciferase reporter assay system that includes, in addition to ERR β Reporter Cells, two optimized media for use during cell culture and in diluting the user's test samples, the reference inverse-agonist 4-Hydroxy Tamoxifen, reagents to prepare Luciferase Detection Reagent, a cell culture-ready assay plate, and a detailed protocol.

Ordering Information

[Order Online »](#)

NGO-IB08011-0096	96 wells
NGO-IB08011-0032	3x32 assays
NGO-IB08012-0384	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 Estrogen-related receptor beta



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