

INDIGlo Luciferase Detection Reagent Kit

INDIGlo Luciferase Detection Reagent (LDR) is formulated to deliver steady, long-lived luminescence signal from mammalian reporter cells expressing firefly luciferase. For convenience, the reagent is provided 'ready to use', without the requirement to first re-constitute powder-form components.

INDIGlo LDR works equally well in quantifying expressed luciferase activity from 96- or 384-well format reporter assays. The resulting light emission is stable over a 2 hr period (Figure 1), thereby making INDIGlo LDR an ideal, economical detection reagent for both low- and high-throughput batch processing of assay plates.

Ordering Information

[Order Online »](#)

NGO-LDR-10	10ml
NGO-LDR-25	25ml
NGO-LDR-50	50ml

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Contents INDIGlo LDR, formulated as a 2x-concentrated reagent
Dilution Buffer, for optional use if performing non-homogenous assays



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