

Human:HeLa NHEJ Reporter Cell Line (Includes I-Sce1 Transfection) (125

assays)

This is a HeLa cell based reporter kit to analyze NHEJ in live cells. Specifically this kit will identify agents (drugs, natural products, small molecules, synthetics, miRNAs, and other genes) that affect or impact NHEJ repair. The kit uses GFP as an in situ readout for NHEJ activity in a cell context. The NHEJ process is activated by transfecting with an I-Sce1 expression plasmid. A specific reporter based assay for NHEJ can be very beneficial for anti-cancer drug discovery projects, learning more about the process of NHEJ repair and establishing intersecting pathways, druggable pathway targets, aging and disease modeling in cells.

Ordering Information

[Order Online »](#)

TOP-DR5001-HNHela- Sce3	125 assays
----------------------------	------------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name NHEJ Based HeLa Requires I-Sce1 Transfection (250ug DNA)



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