

Custom Cell Based Screening Kit for NHEJ (25 assays)

Allows Customer to Create a Cell Line Host for a Non-homologous End Joining (NHEJ)

Reporter This kit will allow you to analyze NHEJ DNA repair in a specific cell lineage or cell line using a neutral GFP gene reporter.

Ordering Information

[Order Online »](#)

TOP-DR5000-Custom-Scce	25 assays
------------------------	-----------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Customer creates cell line hosting NHEJ Reporter (I-Sce Transfectn req.) Includes iHNHeLa cell line control (50ug 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