

# Human:HeLa NHEJ Reporter Cell Line (Includes I-Sce1 Transfection) (25

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-Sce	25 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 (50ug DNA)



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)