

# Custom Cell Based Screening Kit for NHEJ (50 assays)

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

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- Sce2	50 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 (100ug 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)