

# Human Neuroblastoma cell HNSH NHEJ Reporter with I-Sce1 DNA (125

assays)

This is a cell-based reporter kit designed to study NHEJ repair in a human neuroblastoma line (SH-SY-5Y) allowing the customer to screen or identify agents (drugs, natural products, small molecules, synthetics, miRNAs, and genes) that affect or impact the process of DNA repair. The kit uses GFP as an in vivo readout for the NHEJ pathway.

**Cell Based Screening Kit for Non-homologous End Joining (NHEJ) Repair Pathway in a Human Neuroblastoma Cell**

**Ordering Information**

[Order Online »](#)

TOP-DR5000-HNSH-Sce3	125 assays
----------------------	------------

**Manuals, SDS & CofA**

[View Online »](#)

## Handling & Storage

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** NHEJ Based in Human Neuroblastoma Line requires I-Sce1 transfection (250ug 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)