

# Mouse: HNCN1.4-Sce NHEJ Reporter Cell line (Includes I-Sce1 Transfection) (125

assays)

This is a cell-based reporter kit designed to study NHEJ repair in a mouse neuronal cell line that can be differentiated, in vitro. This facilitates identifying 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.  
**Mouse Neuronal Cell Based NHEJ Reporter  
(Requires I-Sce1 Transfection)**

## Ordering Information

[Order Online »](#)

TOP-DR5000-HNCN1-4- Sce3	125 assays
-----------------------------	------------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** NHEJ Based in Mouse Neuronal CN1.4 cells 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)