

Mouse: HNCN1.4-Sce NHEJ Reporter Cell line (Includes I-Sce1 Transfection) (25

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- Sce	25 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 (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