

Mouse: Neuronal Lineage HR Reporter Cell line (Includes I- Sce1 Transfection) (50 assays)

This is a cell-based reporter kit designed to study homologous recombination or HR repair in a mouse neuronal cell line (CN1.4). The cell line expresses numerous neuronal markers, including neurofilaments, glutamic acid decarboxylase 67, neuron-specific enolase, and the BG21 isoform of the myelin basic protein gene. These cells host a reporter gene (GFP) as an in vivo readout for the HR pathway. The cells can be terminally differentiated in vitro.

Ordering Information

[Order Online »](#)

TOP-DR3000-HRCN1-4- Sce2	50 assays
-----------------------------	-----------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name HR based Kit in Mouse Neuronal Cells Requiring I-Sce1 Transfection (100ug 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