

Mouse:3T3 Fibroblast NHEJ Reporter (Includes I-Sce1 Transfection) (125

assays)

DNA Repair pathways in animal cells can be divided into two main categories: HR and NHEJ. HR or homologous recombination is a minor pathway but very important in protecting cells from genotoxicity. NHEJ is the major repair pathway in all animal cells and operates continuously during the entire cell cycle. NHEJ is both error prone but also pivotal in survival of cells since it is a re-ligation pathway that maintains genome integrity.

Ordering Information

[Order Online »](#)

TOP-DR5000-HN3T3-Sce3	125 assays
-----------------------	------------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name NHEJ Based Mouse 3T3 cells Requires I-Sce1 Transfection (250ug 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