

# Mouse:3T3 Fibroblast NHEJ Reporter (Includes I-Sce1 Transfection) (50

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-Sce2	50 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 (100ug 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)