

# Human: Neuroblastoma HR Reporter Cell line (SH- SY5Y, Includes I-Sce1

## Transfection) (125 assays)

This is a cell-based DNA Repair HR reporter that is hosted in human neuroblastoma line SH-SY5Y (or SH cells). The kit is designed to allow the customer to screen or identify agents (drugs, natural products, small molecules, synthetics, miRNAs, pathway, signaling or response genes) that affect or impact the process of HR. The kit uses GFP as an in vivo readout for the HR pathway in a neuro-derived line that can be differentiated. HR is initiated by transfection with an I-Sce1 Expression plasmid (included). Analysis of pooled transfected cells reveals GFP+ cells in 2-3d post-transfection. GFP+ cells have successfully completed HR and can be tracked with single cell resolution.

### Ordering Information

[Order Online »](#)

TOP-DR3000-HRSH-Sce3	125 assays
----------------------	------------

### Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** HR in human Neuroblastoma Cells req. 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)