

# Mouse: 3T3 HR Reporter Cell Line (Includes I-Sce1 Transfection) (50

assays)

This is a mouse 3T3 cell-based reporter kit 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. These cells are easily synchronized and provide the opportunity to analyze HR in the cell cycle.

## Ordering Information

[Order Online »](#)

TOP-DR3000-HR3T3-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 3T3 Cells Requiring 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)