

Mouse: 3T3 HR Reporter Cell Line (Includes I-Sce1 Transfection) (125

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-Sce3	125 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 (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