

Human: HeLa Reporter Cell Line (Includes I-Sce1 Transfection) (50

assays)

This is a HeLa 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. 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-HRHeLa- Sce2	50 assays
----------------------------	-----------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name HR based Kit in HeLa Requiring I-Sce1 Transfection (100ug 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