

Custom Cell Based Screening Kit for (HR) Pathway – no cell line included (50 assays)

Custom cell based screening kit for homologous repair (HR) pathway. This kit allows customer to create a cell line host for HR.

A Kit to Stably Engineer the Homologous Repair (HR) Reporter System in Any Cell Line

Ordering Information

[Order Online »](#)

TOP-DR3000A-Custom- Sce2	50 assays
-----------------------------	-----------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Customer makes own cell line hosting HR Reporter (I-Sce Transfectn req.) No HRHeLa cell line control (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