

Homologous Recombination (HR) Repair Reporter Kit (Transient Assay- Based Kit) (1000 µg)

A kit to interrogate Homologous Repair (HR) Reporter System in Animal Cell Lines: Use of Pooled HR Cells with a GFP Reporter

Customized: You Can Interrogate HR
Ordering Information [Order Online »](#)
Process in A Cell Line Of Your Choosing

TOP-DR3050-4	1000µg
--------------	--------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name HR Pooled Cell analysis: 1000ug DR-HRGFP, 10000ug pSce1, 25ug e-GFP, Puromycin



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