

# Homologous Recombination (HR) Repair Reporter Kit (Transient Assay- Based Kit) (50 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-DR3050-1	50µg
--------------	------

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** HR Pooled Cell analysis: 50ug DR-HRGFP, 50ug pSce1, 25ug e-GFP, Puromycin



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)