

# Human: HeLa HR Reporter Cell line (Includes CRISPR-CAS Transfection)

(125 assays)

This is a cell-based reporter kit designed to study homologous recombination or HR repair in a mouse neuronal cell line (CN1.4). The cell line expresses numerous neuronal markers, including neurofilaments, glutamic acid decarboxylase 67, neuron-specific enolase, and the BG21 isoform of the myelin basic protein gene. These cells host a reporter gene (GFP) as an in vivo readout for the HR pathway. The cells can be terminally differentiated in vitro.

**Cell Based Screening Kit Homologous Repair (HR) Pathway in a Neuroblastoma Cell Line (SH-SY5Y)**

## Ordering Information

[Order Online »](#)

TOP-DR3000-HRHeLa-CPR3	125 assays
------------------------	------------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** HR based Kit in HeLa Requiring CRISPR Transfection (250ug DNA)



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)