

Human P- Glycoprotein / MDR1 Drug Interaction Assay

The materials provided in the assay kit for MDR1 / Human P-Glycoprotein Drug Interaction are formatted to allow for two independent 48-well assay setups. Each aliquot of HCT-Pgp Cells is provided as a single-use assay reagent and provides sufficient volume to dispense 48 assay wells of HCT-Pgp cells into one 96-well plate. If desired, the two aliquots of HCT-Pgp cells may be thawed, combined to generate a single cell suspension, and then dispensed into all wells of an assay plate.

The kit contains two aliquots of HCT-Pgp cells and two black, sterile, collagen-coated 96-well assay plates. In addition to HCT-Pgp cells and assay plates, this kit includes Cell Recovery Medium (CRM-p) used to perform a rapid thaw of the -80°C HCT-Pgp cells, Compound Screening Medium supplemented with daunorubicin (CSM+DR) used to prepare drug treatment media, Wash Buffer, two aliquots of Lysis Reagent, and the potent reference inhibitor Tariquidar.

Ordering Information

[Order Online »](#)

NGO-HPGP-0048	2x48 assays
---------------	-------------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Shipping

Dry Ice

Regulatory Status RUO - Research Use Only

Product Details



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)