

MULTISCREEN™ CHO DHFR- DIVISION ARRESTED CELL LINE

Data sheet

PRODUCT INFORMATION

Catalog Number: DCD0001-1

Lot Number: DCD0001-1-030223

Quantity: 1 vial (4×10^6) frozen cells

Freeze Medium: Cellbanker 2
(Amsbio)

Host cell: CHO dhfr

Recommended Storage: Liquid
nitrogen upon receiving

Propagation Medium: Alpha-MEM,
10% FBS

Application: Functional assays

Figure 1

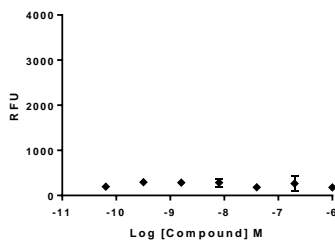


Figure 1: Dose-dependent calcium flux upon treatment with ligand, measured with MULTISCREEN™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01).

FOR RESEARCH USE ONLY.

Multispan Inc. All rights reserved. No part of this document may be reproduced in any form without prior permission in writing.