

MULTISCREEN™ CHO-K1 NFAT DIVISION ARRESTED CELL LINE

Data sheet

PRODUCT INFORMATION

Catalog Number: DCK0001-1

Lot Number: DCK0001-1-032116

Quantity: 1 vial (4 x 10⁶) frozen cells

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: CHO-K1 NFAT

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM/F12, 10% FBS

Stability: Stable for 1-2 days after thawing

Application: Functional assays

Figure 1

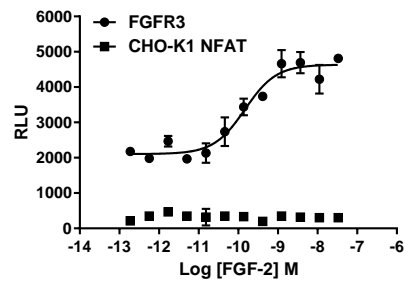


Figure 1: Ligand-induced dose-dependent luciferase activity measured by FlexStation.

FOR RESEARCH USE ONLY.

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