

## MULTISCREEN™ CHO-K1 NFAT CELL LINE

### Data sheet

#### PRODUCT INFORMATION

**Catalog Number:** CK0001-1

**Lot Number:** CK0001-1-032116

**Quantity:** 1 vial ( $4 \times 10^6$ ) frozen cells

**Freeze Medium:** Sigma Freezing Medium (C-6164)

**Host cell:** CHO-K1 NFAT

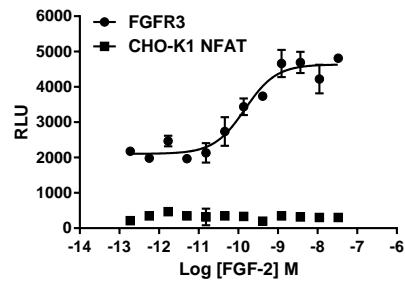
**Recommended Storage:** Liquid nitrogen upon receiving

**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.