

**MULTISCREEN™ DIVISION-ARRESTED CELL LINE  
CHO-K1**

---

**Data sheet**

**PRODUCT INFORMATION**

**Catalog Number:** DC0001-1

**Lot Number:** DC0001-1-072717

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

**Freeze Medium:** Cell banker 2

**Host cell:** CHO-K1

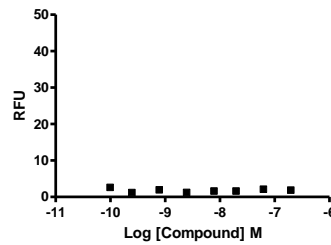
**Recommended Storage:** Liquid nitrogen upon receiving

**Propagation Medium:** DMEM/F12, 10% FBS

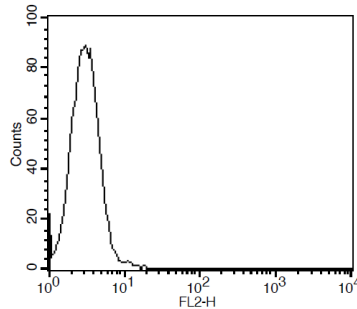
**Stability:** Stable for 1-2 days after thawing

**Application:** Functional assays

**Figure 1**



**Figure 2**



**Figure 1.** Dose-dependent stimulation of calcium flux upon treatment with ligand, monitored with FlexStation. **Figure 2.** Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells.

**FOR RESEARCH USE ONLY.**

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