

**MULTISCREEN™ DIVISION-ARRESTED CELL LINE**  
**CHO-K1 Ga16 GaQI5**

**Data sheet**

**PRODUCT INFORMATION**

**Catalog Number:** DCG0003-1

**Lot Number:** DCG0003-1-053118

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

**Freeze Medium:** Amsbio Cellbanker 2 (11891)

**Host cell:** CHO-K1 Ga16 GaqI5

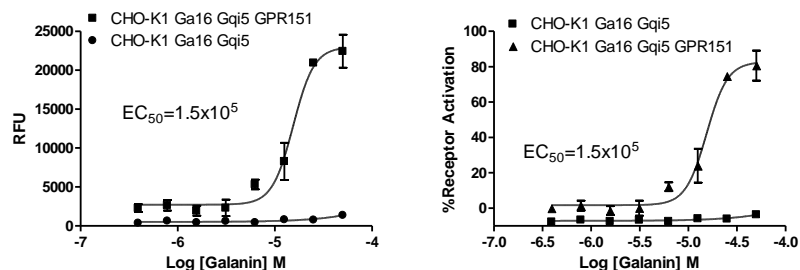
**Recommended Storage:** Liquid nitrogen upon receiving

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

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

**Application:** Functional assays

**Figure 1**



**Figure 1:** Dose-dependent stimulation of calcium flux upon treatment with ligand, monitored with FlexStation.

**FOR RESEARCH USE ONLY.**

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