

MULTISCREEN™ CHO-K1 Gα16 GαQI5 STABLE CELL LINE

Data sheet

PRODUCT INFORMATION

Catalog Number: CG0003-1

Lot Number: CG0003-1-120417

Quantity: 1 vial (2 x 10⁶) frozen cells

Freeze Medium: Cell Banker 2
(Amsbio 11891)

Host cell: CHO-K1 Gα16 GqI5

Recommended Storage: Liquid
nitrogen upon receiving

Propagation Medium: DMEM, 10%
FBS, 250 µg/mL hygromycin, 1 mg/mL
G418

Stability: In Progress

Application: Functional assays

Figure 1

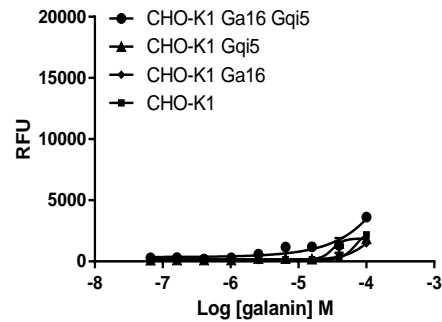


Figure 1. Dose-dependent stimulation of calcium flux upon treatment with ligand, measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01).

FOR RESEARCH USE ONLY.

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