

MULTISCREEN™ CHO-K1 GαQI5 PARENTAL STABLE CELL LINE

Data sheet

PRODUCT INFORMATION

Catalog Number: CG0001-1

Lot Number: CG0001-1-031715

Quantity: 1 vial (2×10^6) frozen cells

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: CHO-K1 Gαq5

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM/F12, 10% FBS

Stability: In progress

Application: Functional assays

Figure 1

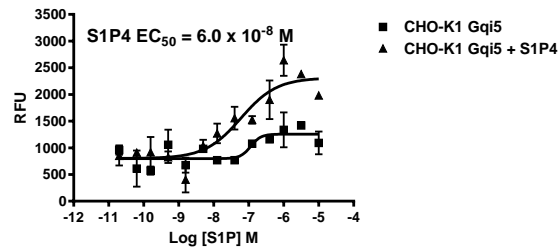


Figure 1: Dose-dependent stimulation of calcium flux upon treatment with ligand, measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01).

FOR RESEARCH USE ONLY.

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