

MULTISCREEN™ DIVISION-ARRESTED CELL LINE CHO-K1 GαQI5

Data sheet

PRODUCT INFORMATION

Catalog Number: DCG0001-1

Lot Number: DCG0001-1-050918

Quantity: 1 vial (4 x 10⁶) frozen cells

Freeze Medium: Cellbanker 2

Host cell: CHO-K1 GαqI5

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM/F12, 10% FBS

Stability: Stability for 1-2 days after thawing

Application: Functional assays

Figure 1

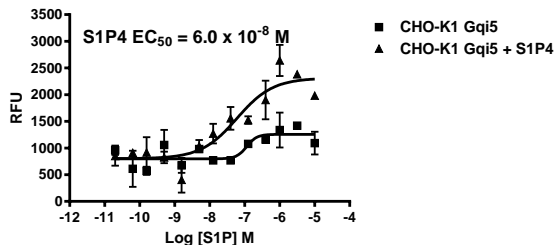


Figure 1: Dose-dependent stimulation of calcium flux upon treatment with ligand, measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01).

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.