

MULTISCREEN™ STABLE CELL LINE HEK293T GαQI5

Data sheet

PRODUCT INFORMATION

Catalog Number: CG0001

Lot Number: CG0001-031915

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

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: HEK293T GαqI5

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM, 10% FBS

Stability: Stable for a minimum of 2 months in continuous culture

Application: Functional assays

Figure 1

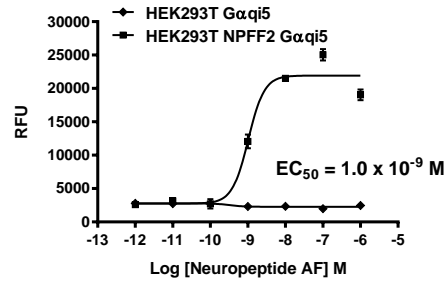


Figure 1: Dose-dependent stimulation of calcium flux upon treatment with ligand, measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01).

FOR RESEARCH USE ONLY.

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