

**MULTISCREEN™ DIVISION-ARRESTED CELL LINE
HUMAN RECOMBINANT APJ RECEPTOR**

PRODUCT INFORMATION

Catalog Number: DC1196a

Lot Number: DC1196a-030615

Quantity: 1 vial (4×10^6) frozen cells

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: HEK293T

Transfection: Expression vector containing full-length human APLNR cDNA (GenBank Accession Number NM_005161) with FLAG tag sequence at N-terminus

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM, 10% FBS

Stability: Stable for 1-2 days after thawing

Data sheet

Background: APJ (also called AGTRL1, angiotensin II receptor-like 1) is a G-protein-coupled receptor that mediates diverse physiological effects of the neuropeptide apelin. It plays a role in the central and peripheral regulation of the cardiovascular system, in water and food intake, and possibly in immune function. APJ is highly expressed in the cardiovascular system and exerts the hypotensive effect *in vivo* and plays a counter-regulatory role against the repressor action of angiotensin II. APJ is also found to be a coreceptor for the entry of several HIV-1 and SIV strains, and apelin blocks the entry of HIV-1 and HIV-2.

Application: Functional assays

Figure 1

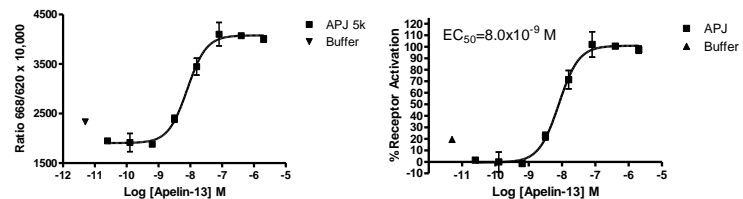


Figure 1. Dose-dependent inhibition of forskolin-stimulated intracellular cAMP level upon treatment with ligand, measured with Multiscreen™ TR-FRET cAMP 1.0 No Wash Assay Kit (Multispan MSCM01).

References:

Cayabyab *et al.* (2000) Apelin, the natural ligand of the orphan seven-transmembrane receptor APJ, inhibits human immunodeficiency virus type 1 entry. *J Virol* 74:11972-11976.

Katugampola *et al.* (2002) Discovery of recently adopted orphan receptors for apelin, urotensin II, and ghrelin identified using novel radioligands and functional role in the human cardiovascular system. *Can J Physiol Pharmacol* 80:369-374.

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.