

**MULTISCREEN™ STABLE CELL LINE**  
**HUMAN RECOMBINANT GPR120 RECEPTOR**

**Data sheet**

**PRODUCT INFORMATION**

**Catalog Number:** CA1522

**Lot Number:** CA1522-061019

**Quantity:** 1 vial (2 x 10<sup>6</sup>) frozen cells

**Freeze Medium:** CellBanker 2 (Amsbio 11891)

**Host cell:** HEK293T

**Transfection:** Expression vector containing full-length human GPR120 cDNA (GenBank Accession Number: NM\_181745.3) with FLAG tag sequence at N-terminus and ARRB2 cDNA (GenBank Accession Number NM\_004313.3)

**Recommended Storage:** Liquid nitrogen upon receiving

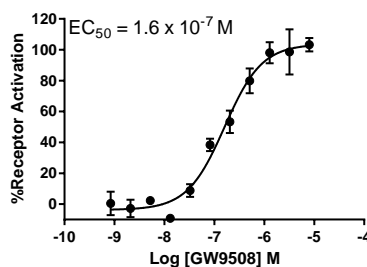
**Propagation Medium:** DMEM, 10% FBS, 1 µg/mL puromycin, 250 µg/mL hygromycin

**Stability:** In Progress

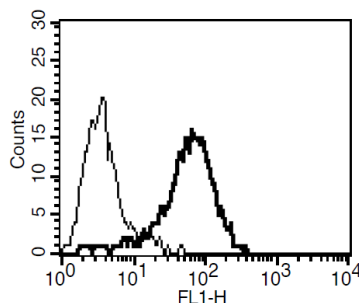
**Background:** GPR120 is a G protein-coupled receptor for the long-chain free fatty acids. GPR120 mediated calcium mobilization, Erk1/Erk2 activation and GLP1 secretion. Unsaturated long-chain FFAs had a dose-dependent stimulatory effect, and  $\alpha$ -linolenic acid was the most potent. GPR120 and GLP1 colocalized in human colonic intraepithelial neuroendocrine cells, and GPR120 may mediate dietary FFA-stimulated GLP1 secretion.

**Application:** Functional assays

**Figure 1**



**Figure 2**



**Figure 1.** Dose-dependent stimulation from arrestin recruitment upon treatment with ligand, monitored on Flexstation III. **Figure 2.** Receptor Expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells; thick line: receptor-expressing cells.

**References:**

Fredriksson *et al.* (2003) Seven evolutionarily conserved human rhodopsin G protein-coupled receptors lacking close relatives. *FEBS Lett* 554:381-388.

Hirasawa *et al.* (2005) Free fatty acids regulate gut incretin glucagon-like peptide-1 secretion through GPR120. *Nature Med* 11:90-94.

**FOR RESEARCH USE ONLY.**

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