

**MULTISCREEN™ STABLE CELL LINE  
HUMAN RECOMBINANT GPR81 RECEPTOR**

**PRODUCT INFORMATION**

**Catalog Number:** C1134

**Lot Number:** C1134-110617

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

**Freeze Medium:** Cell Banker 2  
(Amsbio 11891)

**Host cell:** HEK293T

**Transfection:** Full-length Human GPR81 (GenBank Accession Number NM\_032554.3) with FLAG-tag sequence at the N-terminus

**Recommended Storage:** Liquid nitrogen upon receiving

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

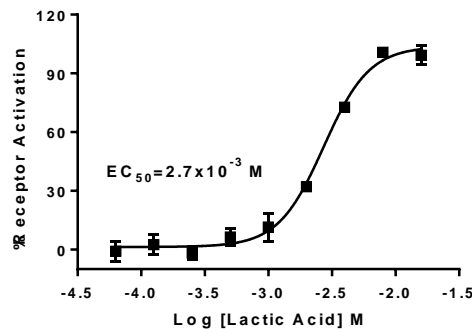
**Stability:** In Progress

**Data sheet**

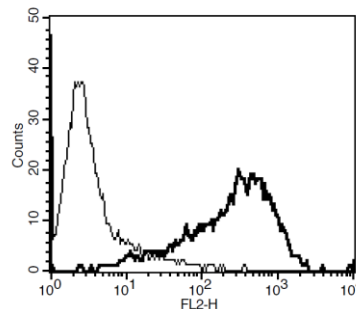
**Background:** GPR81 is an orphan G protein-coupled receptor homologous to GPR109A and GPR109B. It is highly expressed in adipose tissues and also regarded as a sensor for Lactate and can mediate an anti-lipolytic effect of lactate. GPR81 can modulate the free fatty acid pool under exercise or conditions of oxygen deficit and is target for the treatment of dyslipidemia and other metabolic disorders.

**Application:** Functional assays

**Figure 1**



**Figure 2**



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

Ge *et al* (2008) Elucidation of signaling and functional activities of an orphan GPCR, GPR81. *Journal of Lipid Research* 49:797-803.

Liu *et al.* (2009) Lactate Inhibits Lipolysis in Fat Cells through Activation of an Orphan G-protein-coupled Receptor, GPR81. *Journal of Biological Chemistry* 284: 2811–2822.

Cai *et al* (2008) Role of GPR81 in lactate-mediated reduction of adipose lipolysis. *Biochem Biophys Res Commun*, 377(3): 987-991

**FOR RESEARCH USE ONLY.**

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