

**MULTISCREEN™ DIVISION ARRESTED CELL LINE  
HUMAN RECOMBINANT D2 RECEPTOR**

**Data sheet**

**PRODUCT INFORMATION**

**Catalog Number:** DC1336

**Lot Number:** 031411

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

**Freeze Medium:** Sigma Freezing Medium (C-6164)

**Host cell:** HEK293T

**Transfection:** Expression vector containing full-length human DRD2 cDNA (GenBank accession number NM\_000795.3) with FLAG tag sequence at N-terminus

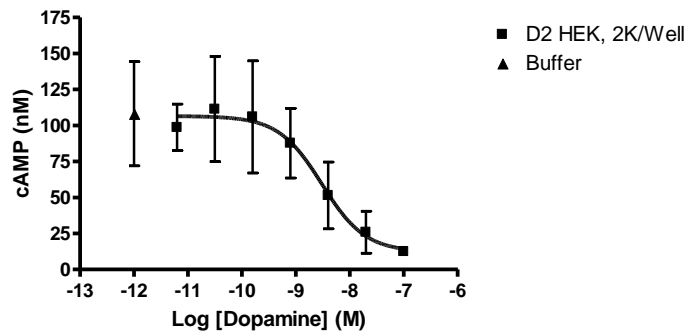
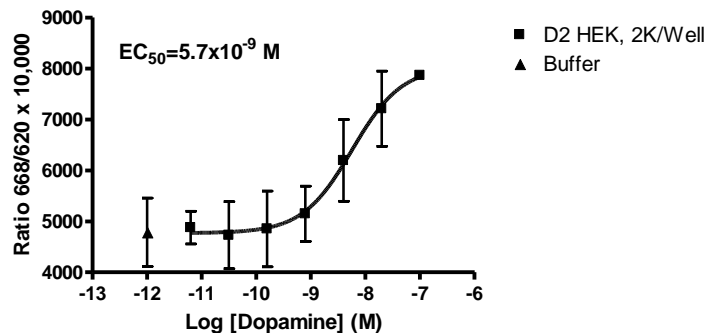
**Recommended Storage:** Liquid nitrogen upon receiving

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

**Stability:** Stable for 1-2 days after thawing

**Background:** The human dopamine receptor DRD2 (D2) is a G protein-coupled receptor for dopamine. It can be found on postsynaptic dopaminergic neurons that are centrally involved in reward-mediating mesocorticolimbic pathways. Signaling through dopamine D2 receptors governs physiological functions related to locomotion, hormone production, and drug abuse.

**Application:** Functional assays



**References:**

Grandy *et al.* (1989) Cloning of the cDNA and gene for a human D2 dopamine receptor. *Proc Nat Acad Sci USA* 86:9762-9766.

Neville *et al.* (2004) Identification and characterization of ANKK1: a novel kinase gene closely linked to DRD2 on chromosome band 11q23.1. *Hum Mutat* 23:540-545.

Usiello *et al.* (2000) Distinct functions of the two isoforms of dopamine D2 receptors. *Nature* 408:199-203.

**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.