



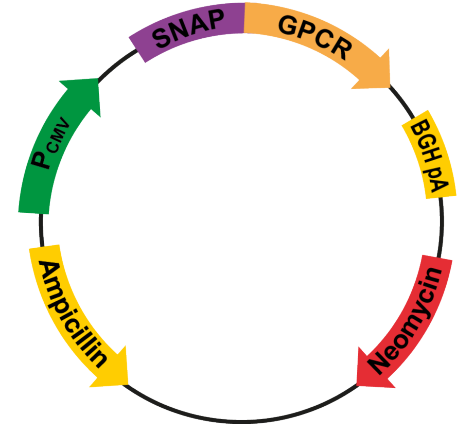
# Dopamine PLASMID TAG-LITE® pSNAP-D2

## Receptor: Dopamine receptor type 2 isoform short (Variant 2)

PRODUCT INFORMATION - rev 1 (Feb. 2010)

**Part#:** PSNAPD2  
**Storage temperature:** -20°C  
**Quantity:** 10 µg/ 10 µL H2O

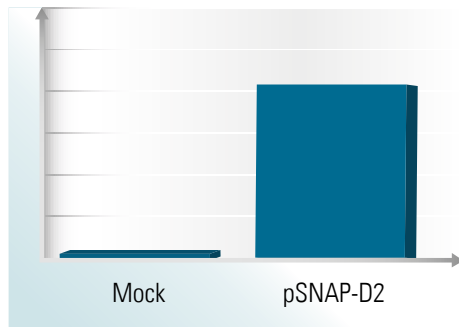
**Plasmid name:** pSNAP-D2  
**Plasmid size:** 7283 bp  
**D2-ORF size:** 1914 bp  
**ORF sequence\*:** NM\_016574.2  
**\*based on NCBI nucleic sequence accession number**  
**Promoter:** CMV  
**Host:** Mammalian  
**Antibiotic resistance for E. Coli:** Ampicillin  
**Mammalian selection marker:** Neomycin



Tag-lite pSNAP-D2

For research use only. Not for use in diagnostic procedures.

### QUALITY CONTROL



Plasmid	Normalized 620 nm signal*
pSNAP-D2	24

\*The plasmid is transiently expressed into HEK 293. Cells are plated and labelled 24 hours after transfection with the Tag-lite SNAP-Lumi4™-Tb (SSNPTBC) according to the protocol described on the web site. Signal is measured at 620 nm and the value normalized versus the value obtained with a mock plasmid expressed as negative control.

### SEQUENCE & MSDS

Files can be downloaded from Tag-lite section on the website [www.cisbio.com](http://www.cisbio.com).

Sequences are available under a txt format

#### FOR MORE INFORMATION

Europe and other countries +33(0)466 796 705 U.S. and Canada 1-888-963-4567

China +86-21-5018-9880 Japan +81-(0)43-306-8712

Visit [www.cisbio.com](http://www.cisbio.com) to find a list of our regional distributors