## Dopamine receptor D<sub>2</sub> polyclonal antibody

Dopamine is the predominant catecholamine neurotransmitter in the mammalian brain, where it regulates locomotor, emotional, memory, and dietary intake functions, as well as peripheral roles such as hormone secretion and renal function. Dopamine receptors (D1-D5) are members of the seven transmembrane domain containing GPCR family, and are the target of several anti-psychosis and anti-Parkinsonian drugs. The D2, D3, and D4 receptors signal via coupling of Gi/Go alpha G-protein subunits to inhibit adenylyl cyclase activity and modulate cellular potassium and calcium ion channel conductance. Expression of the D2 receptor is found primarily in the brain (striatum, olfactory tubercle, and nucleus accumbens) and in the anterior and intermediate lobes of the pituitary.

This antibody is covered by our Worry-Free Guarantee.

Citations: 4

View Online »

**Ordering Information** 

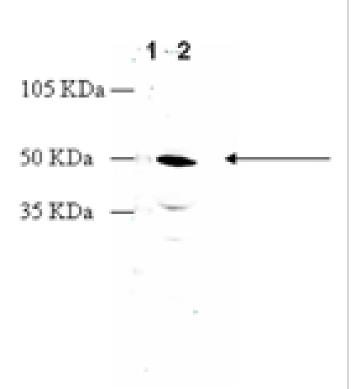
**Order Online** »

ADI-905-740-100

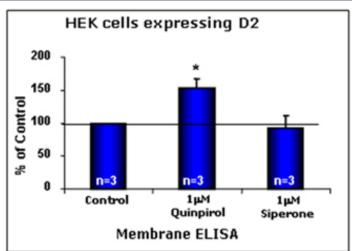
100µg

Manuals, SDS & CofA

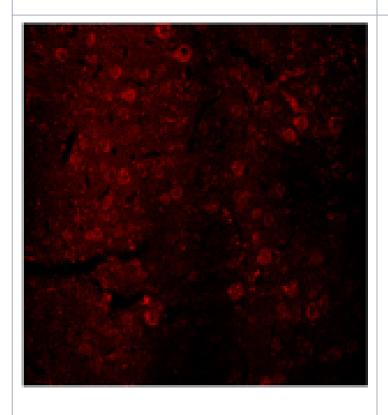
**View Online »** 



Western blot analysis: MW marker (1) and 10µg of rat heart membrane extract (2) probed with Dopamine Receptor (D2) pAb at 4.0µg/ml.



Cellular ELISA: HEK293 cells expressing dopamine receptor (D2) were treated with indicated concentrations of agonist (Quinpirol) and antagonist (Siperone) and probed with Dopamine Receptor (D2) pAb (1:1000 of a stock solution of  $1\mu g/\mu L$ ) by cellular ELISA. Data from vehicle treated cells were taken as 100%. Results are the mean  $\pm$ SEM (n=3).



Immunohistochemistry analysis: Fluorescent confocal photomicrograph of rat cerebral cortex. Neurons are labeled with Dopamine Receptor (D2) pAb in conjunction with Cy3 conjugated rabbit IgG antibody (H+L).

## **Handling & Storage**

**Handling** Keep on ice at all times.

Long Term Storage -20°C

Shipping Blue Ice

## Regulatory Status RUO - Research Use Only

## **Product Details**

Alternative Name DRD, D2R

**Application** ELISA, IHC, WB

**Application Notes**Suitable for cell-based and membrane-based ELISA.

Predicted MW of ~51kDa.

Crossreactivity Predicted species reactivity based on epitope sequence

homology: rhesus macaque, chimpanzee, horse, arctic

fox, dog, opossum, ferret.

Formulation Liquid. In PBS containing 50% glycerol and 0.01% sodium

azide.

**Host** Rabbit

**Immunogen** Synthetic peptide corresponding to the sequence near the

N-terminus of mouse dopamine receptor D2</sub.

Purity Detail Peptide affinity purified.

Species Reactivity Human, Mouse, Rat

UniProt ID P61168

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

