

CP 55,940

CP 55,940 is a synthetic cannabinoid molecule developed by Pfizer that has similar but more potent effects to THC in behavioral and receptor binding tests. It is a cannabinoid receptor agonist, with EC₅₀=0.2, 0.3, and 5nM for CB1, CB2, and GRP55 respectively. Currently used to study the endocannabinoid system.

Materials Provided

Catalog Number:	92-1343
Quantity Supplied:	1 vial(s), 5 mg per vial

Description

DiscoverX control ligands have been functionally tested and validated for optimal use with all cell line targets and Assay Ready kits.

Product Information

Molecular Weight:	376.6
Source:	Synthetic
Purity:	>98%
Endotoxin Level:	N/A
Formulation:	N/A
Storage:	-20°C Please avoid multiple freeze/thaw cycles.

Reagent Preparation

To avoid condensation, equilibrate the vial to ambient temperature before opening.

Stock Concentration:	10 mM
Reconstitution Volume:	1.330 ml
Reconstitution Solvent:	DMSO

Handling Instructions:

Soluble in ethanol, DMSO and DMF all at ~30 mg/ml. Sparingly soluble in aqueous buffers, ethanol:PBS (1:3, pH 7.2, 0.25 mg/ml).

For Research Use Only

Ordering: +1.510.979.1415 option 4 or e-mail CustomerServiceDRX@eurofins.com
Technical support: +1.510.979.1415 option 5 or e-mail DRX_SupportUS@eurofinsUS.com
General product information: www.discoverx.com