

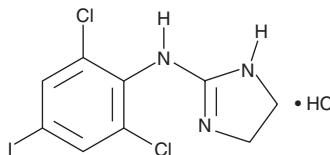
PRODUCT INFORMATION



p-iodo-Clonidine (hydrochloride)

Item No. 17177

CAS Registry No.: 108294-57-1
Formal Name: N-(2,6-dichloro-4-iodophenyl)-4,5-dihydro-1H-imidazol-2-amine, monohydrochloride
MF: C₉H₈Cl₂IN₃ • HCl
FW: 392.5
Purity: ≥98%
UV/Vis.: λ_{max}: 216, 247 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

p-iodo-Clonidine (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the *p*-iodo-clonidine (hydrochloride) in the solvent of choice, which should be purged with an inert gas. *p*-iodo-Clonidine (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of *p*-iodo-clonidine (hydrochloride) in ethanol and DMSO is approximately 1 mg/ml and approximately 0.25 mg/ml in DMF.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of *p*-iodo-clonidine (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of *p*-iodo-clonidine (hydrochloride) in PBS (pH 7.2) is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

p-iodo-Clonidine is a partial agonist of the α₂-adrenergic receptor, minimally inhibiting adenylate cyclase through this Gi protein-coupled receptor although it binds avidly (K_i = 1.0 nM).¹ It potentiates platelet aggregation by ADP (EC₅₀ = 1.5 μM) and inhibits epinephrine-induced aggregation (IC₅₀ = 5.1 μM).¹

Reference

1. Gerhardt, M.A., Wade, S.M., and Neubig, R.R. *p*-[125I]iodoclonidine is a partial agonist at the alpha 2-adrenergic receptor. *Mol. Pharmacol.* **38**(2), 214-221 (1990).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/14/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM