Diphenyleneiodonium chloride

NADPH oxidase inhibitor

Binds strongly to flavoproteins and is thus a powerful and specific inhibitor of several important enzymes, including nitric oxide synthase (NOS), NADPH-ubiquinone oxidoreductase, NADPH oxidases and NADPH cytochrome P450 oxidoreductase. Nitric oxide synthase, which shows significant homology with cytochrome P450 reductase, has been shown to be irreversibly inhibited by this compound in cultured peritoneal macrophages (IC_{50} =30nM).

Citations: 24

View Online »

Ordering Information

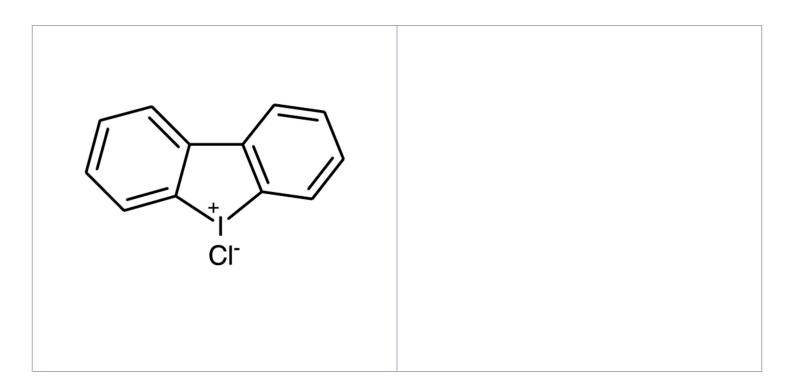
Order Online »

BML-CN240-0010

10mg

Manuals, SDS & CofA

View Online »



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Handling Protect from light.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name DPI chloride, Dibenziodolium chloride

Appearance White to off-white solid.

CAS 4673-26-1

Couple Target Cytochrome P450 reductase, NOS

Couple Type Inhibitor

Formula $C_{12}H_8I$. CI

Identity Determined by 1H-NMR and MS.

MW 314.5

Purity ≥98% (HPLC)

Soluble in DMSO (warm) (5mg/ml).

Last modified: May 29, 2024

