Argatroban

Thrombin inhibitor

Argatroban is a synthetic peptidomimetic thrombin inhibitor which is in clinical use. It presents significant advantages as an anticoagulant in deep vein thrombosis, stroke prevention in atrial fibrillation and thromboembolic disease and especially in patients with coronary artery disease and heparin-induced thrombocytopenia.

Citations: 4

View Online »

Ordering Information

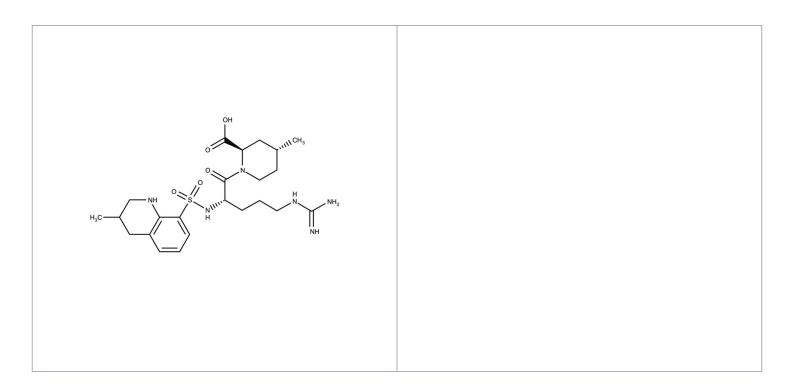
Order Online »

BML-PI146-0010

10mg

Manuals, SDS & CofA

View Online »



Handling & Storage

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

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance White to off-white powder.

CAS 74863-84-6

Couple Target Thrombin

Couple Type Inhibitor

Formula $C_{23}H_{36}N_6O_5S$

Identity Determined by NMR and LCMS.

MW 508.6

Purity ≥98% (HPLC)

Solubility Soluble in DMSO (25mg/ml) or 100% ethanol (25mg/ml).

Last modified: May 29, 2024

