Decadienal

Induces mitochondrial dysfunction and oxidative stress

Decadienal is a lipid peroxidation product which is an aldehydic bifuntional electrophile. It induces mitochondrial dysfunction and oxidative stress possibly via covalent modification of cytochrome c. It can induce cell proliferation via the p27 pathway in human bronchial epithelial cells.

Ordering Information

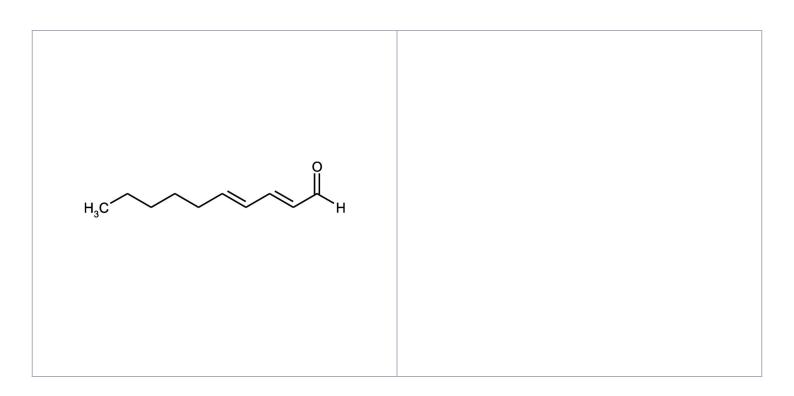
Order Online »

BML-HA005-0020

20mg

Manuals, SDS & CofA

View Online »



Handling & Storage

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

at -20°C for up to 1 year. Store solutions at -20°C for up to 3 months.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Appearance Pale yellow oil.

CAS 25152-84-5

Formula $C_{10}H_{16}O$

MW 152.2

Purity ≥95%

Solubility Soluble at 25mg/ml in DMSO, ethanol, or DMF.

Last modified: May 29, 2024

