

DAF-FMCatalog number: 16296
Unit size: 1 mg**Product Details**

Storage Conditions	Freeze (<-15 °C), Minimize light exposure
Expiration Date	12 months upon receiving

Chemical Properties

Appearance	Solid
Molecular Weight	412.35
Soluble In	DMSO

Applications

DAF-2 is the original compound and is widely used for detecting nitric oxide (NO) in cells. DAF-FM is an improved version of DAF. Compared with DAF-2, DAF-FM is a much more sensitive and photo-stable fluorescent probe for detecting NO. In addition, DAF-FM is much less pH sensitive than DAF-2, essentially pH-independent in the physiological pH range. DAF-FM is not quite as cell-permeable as the diacetate derivative, DAF-FM, but it does not need to be activated by cytosolic enzymes.