

Texas Red–dextran conjugate (average MW = ~20K, fixable)

Catalog Number: 21722

Unit Size: 25 mg

Product Details

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
--------------------	--

Expiration Date	12 months upon receiving
-----------------	--------------------------

Chemical Properties

Appearance	Solid
------------	-------

Molecular Weight	~20,000
------------------	---------

Soluble In	Water
------------	-------

Spectral Properties

Excitation Wavelength	586 nm
-----------------------	--------

Emission Wavelength	603 nm
---------------------	--------

Applications

Texas Red–dextran conjugate (average MW = ~20K, fixable) is a lysine-derivatized dextran intended for fluorescence-based tracing and transport studies in cells, tissues, and whole organisms. Conjugated to Texas Red (Ex/Em = 586/603 nm), it produces a bright, photostable red-orange signal compatible with standard fluorescence microscopy as well as advanced imaging modalities. The 20-kDa dextran backbone provides defined diffusional properties that support applications such as anterograde and retrograde neuronal labeling, endothelial barrier assessment, and measurements of extracellular or intracellular transport.

The presence of primary amine groups allows the conjugate to become covalently crosslinked to proteins during aldehyde fixation, ensuring that the fluorescent signal remains associated with the labeled structures through permeabilization, immunostaining, tissue clearing, and long-term storage. This fixability permits reliable correlation between the tracer distribution and additional molecular or structural markers introduced after fixation.