



## Aminoallyl-UTP-PEG<sub>5</sub>-IR680LT

UTP-PEG<sub>5</sub>-IR680LT

5-(3-Aminoallyl-PEG<sub>5</sub>)-uridine-5'-triphosphate, labeled with IR680LT, Triethylammonium salt

Cat. No.	Amount
NU-821-PEG5-IR680LT	10 µl (5 mM)

**For general laboratory use.**

**Shipping:** shipped on gel packs

**Storage Conditions:** store at -20 °C

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

**Shelf Life:** 12 months after date of delivery

**Molecular Formula:** C<sub>75</sub>H<sub>97</sub>N<sub>6</sub>O<sub>40</sub>P<sub>3</sub>S<sub>6</sub>

**Molecular Weight:** 2007.88 g/mol (free acid)

**Exact Mass:** 2006.33 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

**Form:** solution in 10 mM Tris-HCl

**Color:** blue

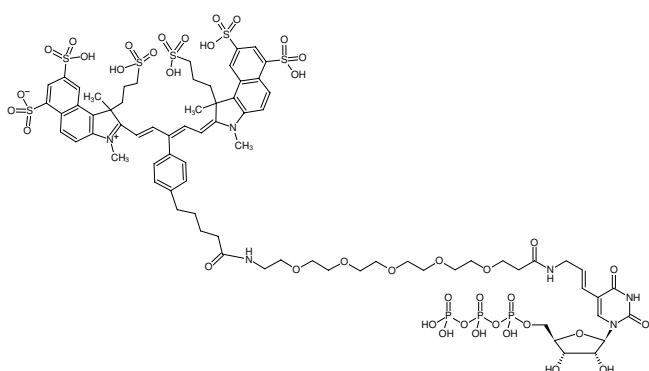
**Concentration:** 5.0 mM - 5.5 mM

**pH:** 7.5 ± 0.5

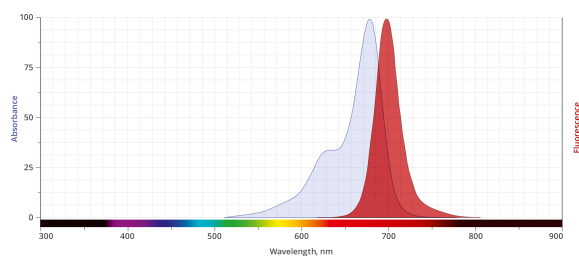
**Spectroscopic Properties:** λ<sub>exc</sub> 679 nm, λ<sub>em</sub> 694 nm, ε 190.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)

**Applications:**

RNA Labeling by in vitro Transcription with T7 RNA Polymerase (in-house tested)



Structural formula of Aminoallyl-UTP-PEG<sub>5</sub>-IR680LT



excitation and emission spectrum of IR680LT