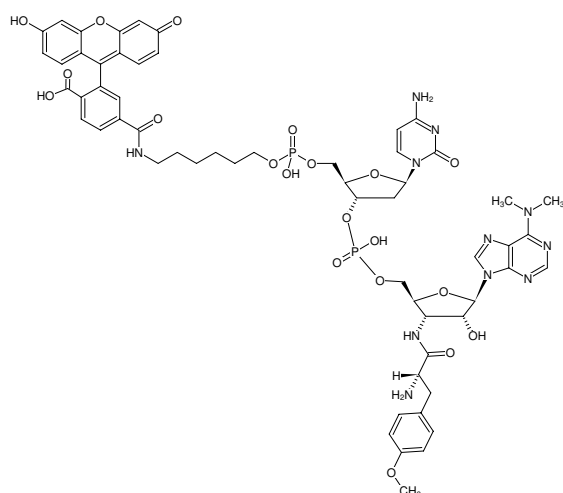


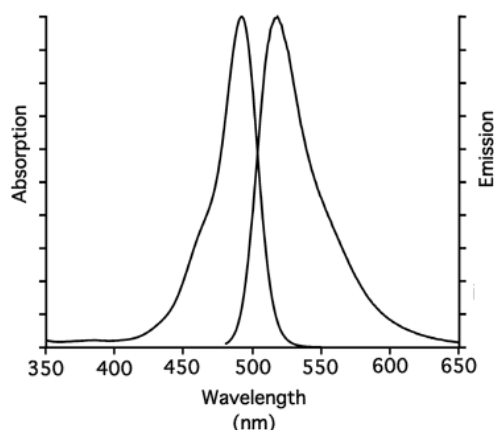


## 6-FAM-dC-puromycin

Cat. No.	Amount
NU-925-6FM-S	100 µl (0,1 mM)
NU-925-6FM-L	5 x 100 µl (0,1 mM)



Structural formula of 6-FAM-dC-puromycin



excitation and emission spectrum of 6-FAM

**For research use only!**

**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>58</sub>H<sub>65</sub>N<sub>11</sub>O<sub>20</sub>P<sub>2</sub> (free acid)

**Molecular Weight:** 1298.15 g/mol (free acid)

**Exact Mass:** 1297.39 g/mol (free acid)

**CAS#:** 436083-76-0

**Purity:** ≥ 95 % (HPLC)

**Form:** solution in water

**Color:** yellow

**Concentration:** 0.10 mM - 0.11 mM

**pH:** 7.5 ± 0.5

**Spectroscopic Properties:** λ<sub>exc</sub> 499 nm, λ<sub>em</sub> 517 nm, ε 83.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 9.0)

**Selected References:**

Starck *et al.* (2004) A general approach to detect protein expression in vivo using fluorescent puromycin conjugates. *Chem. Biol.* **11** (7):999.

Kawahashi *et al.* (2007) High-throughput fluorescence labelling of full-length cDNA products based on a reconstituted translation system. *J. Biochem.* **141** (1):19.