

**AF488 Streptavidin Conjugate**

Streptavidin, Alexa Fluor® 488 Conjugate

Cat. No.	Amount
RNT-015	1 mg

For research use only!**Shipping:** shipped on gel packs**Storage Conditions:** store at -20 °C**Additional Storage Conditions:** store dark**Shelf Life:** 12 months after date of delivery**Form:** solid containing lyophilised PBS buffer (pH 7.2) and 0.1% BSA (without sodium azide)**Color:** yellow to orange**Solubility:** water**Spectroscopic Properties:** λ_{exc} 499 nm, λ_{em} 520 nm, ϵ 73.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)**Description:**

AF488 Streptavidin Conjugate (also known as Alexa Fluor® 488 Streptavidin Conjugate) is ideally suited for the specific, fluorescence-based detection of biotinylated biomolecules (e.g. RNA or DNA) in histochemical, flow cytometry, blotting and microplate assays. The detection of Biotin relies on its high affinity ($K_D = 10^{-15}$ M) to Streptavidin that is labeled with the photostable, hydrophilic AF488 dye (also known as Alexa Fluor® 488). AF488 can be detected with the standard FITC filter set.

Degree of Labeling (molar dye/protein ratio): 2 - 4**Preparation of 1 mg/ml working solution:**Add 1 ml ddH₂O, vortex and spin-down briefly.**Selected References:**

[1] Diamandis *et al.* (1991) The biotin-(strept)avidin system: principles and applications in biotechnology. *Clin. Chem.* **37**:625.