

Conjugation Protocol for Amine Reactive CoraFluor™ Reagents

In Brief

Amine reactive CoraFluor™ reagents contain pentafluorophenyl esters (PFP esters) which can be conjugated to (non-protonated) aliphatic amine groups. The primary reactive species for protein amine-conjugation are the ϵ -amino groups of lysine residues.

To avoid protonating these groups, it is important to perform the reaction at a slightly basic pH. In addition, buffers containing primary amines should be avoided, since they will compete for conjugation with the PFP ester.

Please note that PFP esters can be moisture sensitive, so handle accordingly. Where possible, handle and store CoraFluor™ reagents in the dark.

Conjugation Protocol

1. Prepare a 100 μ L aliquot of an antibody, protein or nanobody at a concentration of ≥ 1 mg/mL in reaction buffer (100 mM sodium bicarbonate buffer, pH 8.5) using a 0.5 mL, 7 kDa molecular weight cutoff (MWCO) Zeba™ spin desalting column (Thermo Fisher 89882) according to the manufacturer's protocol.
2. Add the antibody/protein/nanobody solution to the amine reactive CoraFluor™ to achieve a molar ratio of ~ 5 – $15\times$ CoraFluor™ to antibody/protein, or 4 – $5\times$ CoraFluor™ to nanobody.

If performing multiple conjugations, reconstitute CoraFluor™ in 2.5 mM dry DMSO or DMAc and perform conjugation reactions with a final DMSO or DMAc content $< 10\%$.

The molar equivalents of CoraFluor™ can be adjusted accordingly depending on size of the protein and desired degree of labeling.

3. Briefly vortex the reaction mixture and incubate at room temperature for 1 h.
4. Remove organic solvent and unreacted PFP ester complex by buffer exchange into desired storage buffer (e.g. 50 mM sodium phosphate buffer, pH 7.4, with 150 mM NaCl and 0.05% (vol/vol) TWEEN®-20) using a 0.5 mL, 7 kDa MWCO Zeba™ spin desalting column, according to the manufacturer's protocol.
5. Determine concentration and degree of labeling (DOL) using the below calculations.

Degree of Labeling Calculation

- Determine the corrected absorbance at 280 nm value ($A_{280,corr}$) of the antibody/nanobody/protein conjugate by measuring A_{280} and A_{340} using:

$$A_{280,corr} = A_{280} - (A_{340} \times c.f.)$$

c.f. is the correction factor for the terbium complex contribution to A_{280} and is equal to 0.157.

- Determine the concentration of antibody/protein/nanobody conjugate, c_{ab} (in M) using:

$$c_{ab} = \frac{A_{280,corr}}{\epsilon_{ab}} \times b$$

where ϵ_{ab} is the antibody/protein/nanobody extinction coefficient at A_{280} and b is the path length in centimeters.

- Determine the concentration of covalently bound terbium complex, c_{Tb} (in M) using:

$$c_{Tb} = \frac{A_{340}}{\epsilon_{Tb}} \times b$$

where ϵ_{Tb} is the complex extinction coefficient at A_{340} equal to $22,000 M^{-1} cm^{-1}$, and b is the path length in centimeters.

- Calculate the degree of labeling (DOL) using:

$$DOL = \frac{c_{Tb}}{c_{ab}}$$

TWEEN is a registered trademark of Croda International PLC.

Zeba is a registered trademark of Thermo Fisher Scientific.

References

Payne *et al* (2021) Bright and stable luminescent probes for target engagement profiling in live cells. *Nat.Chem.Biol.* **17** 1168 PMID:34675420.

For research use or manufacturing purposes only. Trademarks and registered trademarks are the property of their respective owners. 7615723165_102024

biotechne // Global Developer, Manufacturer, and Supplier of High-Quality Reagents, Analytical Instruments, and Precision Diagnostics.

INCLUDES R&D Systems™ Novus Biologicals™ Tocris Bioscience™ ProteinSimple™ ACD™ ExosomeDx™ Asuragen® Lunaphore™