
Important Information: Custom Antibody & Protein Conjugates

Unless ordered otherwise, all custom conjugates are delivered as a single aliquot in a 4°C shipment.

Storage and Stability

- Conjugate stability is dependent on the biomolecule and needs to be determined empirically.
 - Typical antibody conjugates are stable for > 1 year at 4°C.
 - Most fluorescent conjugates are stable for at least 6 months when stored properly.
 - Freeze/thaw cycles should be avoided; if -20°C is required, the conjugate solution should be aliquoted and snap frozen to preserve activity.
- Upon receipt, aliquot if desired and store at 4°C for intermediate length storage.
- DO NOT freeze R-Phycoerythrin or Qdot™ conjugates.
- Conjugate expiration is not assigned as stability is dependent on the biomolecule.

Handling Guidelines

- Wear gloves when handling protein conjugates to avoid contamination (enzymatic, bacterial, etc.).
- Centrifuge conjugate before first use. If precipitation occurs, centrifuge conjugate and pipette soluble conjugate from the supernatant.
- Conjugate analytical data is on the provided Certificate of Analysis.
- Functional QC tests are not performed due to protein-specific testing needs. All measures are taken during conjugation to preserve protein activity.
- Protect fluorescent conjugates from direct light:
 - Polymer and tandem dye conjugates are highly sensitive to light exposure.
 - Other fluorescent dye conjugates are mildly to moderately sensitive to light exposure.
- DO NOT add sodium azide to HRP-containing conjugates.

For more information visit: <https://www.thermofisher.com/us/en/home/life-science/antibodies/antibody-labeling/custom-fluorescent-antibody-protein-labeling.html>

Please contact our skilled team of Custom Conjugation Services Project Managers with any questions: lsg-eug-pmoteam@thermofisher.com.