

Protein Standards for Ion Exchange HPLC

Instructions for Reconstitution and Use

Protein Anion Exchange Standard

Catalog #125-0561

Add 2.0 ml of the low salt buffer to be used (refer to operating instructions for specific column). Swirl the vial gently. The lyophilized proteins should dissolve completely. If they don't, filter or centrifuge the standard before using it. The proteins are (in typical order of elution):

1. Myoglobin
2. Conalbumin
3. Ovalbumin
4. Soybean trypsin inhibitor

Unused standard can be stored at 4°C for several days or frozen for up to 1 year from the time of reconstitution.

Protein Cation Exchange Standard

Catalog #125-0562

Add 2.0 ml of the low salt buffer to be used (refer to operating instructions for specific column). Swirl the vial gently. The lyophilized proteins should dissolve completely. If they don't, filter or centrifuge the standard before using it. The proteins are (in typical order of elution):

1. Myoglobin
2. Ribonuclease A
3. Cytochrome c

Unused standard can be stored at 4°C for several days or frozen for up to 1 year from the time of reconstitution.