

Product Information

Fluorescent Molecular Weight Marker (M.W. 20,000–200,000)

Catalog Number **F3526**

Storage Temperature –20 °C

TECHNICAL BULLETIN

Product Description

The Fluorescent Molecular Weight Marker is prepared by conjugation of a highly fluorescent dye to six proteins having a wide range of molecular masses (see Table 1). The conjugates produce sharp bands in SDS-polyacrylamide gel electrophoresis and after transfer to membranes. The Fluorescent Molecular Weight Marker is easily visualized with UV light. The marker may also be visualized with standard Brilliant Blue stains, with quenching of the fluorescence.

Table 1.
Protein Mixture in F3526

Protein	Approximate Molecular Mass (Daltons)
Trypsin Inhibitor, soybean	20,000
Carbonic Anhydrase, bovine erythrocyte	29,000
Alcohol Dehydrogenase, equine liver	39,800
Albumin, bovine serum	66,000
β-Galactosidase, <i>E. coli</i>	116,000
Myosin, rabbit muscle	200,000

Precautions and Disclaimer

This product is for R&D use only, not for drug, household, or other uses. Please consult the Material Safety Data Sheet for information regarding hazards and safe handling practices.

Preparation Instructions

Reconstitute the contents of each vial with 250 µl of water and vortex for a few seconds until completely dissolved. Aliquot and freeze immediately. Store in a container protected from light. Repeated freezing and thawing is **not** recommended.

The resulting protein solution contains 62 mM Tris-HCl, pH 8.0, 1 mM EDTA, 3% sucrose, 0.5% dithiothreitol, 2% SDS, and 0.005% bromophenol blue.

Storage/Stability

Store the product in the dark at –20 °C. After reconstitution, aliquot and freeze. Store in a container protected from light. Repeated freezing and thawing is **not** recommended.

Procedure

1. Incubate the markers in a 65 °C water bath for 5 minutes immediately before loading onto the gel.
Note: Avoid boiling the sample because it may promote degradation of the marker.

2. Recommended sample volumes:

Standard size gel (16 × 14 cm) - 10–20 µl/well

Mini-gel (10 × 10 cm) - 5–10 µl/well

3. During or after electrophoresis, protein bands can be visualized on gels or transfer membranes using a UV lamp or a UV transilluminator.
Note: Due to the large size of myosin, it will not transfer completely to membranes. Band intensity will vary with transfer conditions.

KR,MAM 07/09-1

Sigma brand products are sold through Sigma-Aldrich, Inc.

Sigma-Aldrich, Inc. warrants that its products conform to the information contained in this and other Sigma-Aldrich publications. Purchaser must determine the suitability of the product(s) for their particular use. Additional terms and conditions may apply. Please see reverse side of the invoice or packing slip.