GloMelt™ Thermal Shift Protein Stability Kit

Excitation/Emission

Product attributes 468/507 nm

Call us: 800-304-5357 Email: btinfo@biotium.com

Kit for fluorescence-based measurement of protein unfolding or thermal stability in thermal shift assay, also called Protein Thermal ShiftTM, differential scanning fluorimetry, or Thermofluor assay.

Product Description

The GloMelt™ Thermal Shift Protein Stability Kit is used to detect protein unfolding or measure thermal stability by performing a thermal shift assay, also called Protein Thermal Shift™, differential scanning fluorimetry, or Thermofluor assay.

Also view our GloMelt™ 2.0 Thermal Shift Protein Stability Kit which features a Recombinant Monoclonal Mouse IgG1 for performing control reactions. Compared to the polyclonal goat IgG control provided in the original GloMelt™ kit, the recombinant mouse antibody included in the GloMelt™ 2.0 Kit generates higher signal and a sharper melt peak between for robust assay validation.

About GloMelt™ Thermal Shift Protein Stability Kit

GloMelt™ dye undergoes fluorescence enhancement upon binding to hydrophobic regions of denatured proteins, and therefore can be used to detect protein unfolding or measure thermal stability by performing a thermal shift assay, also called Protein Thermal Shift™, differential scanning fluorimetry, or Thermofluor

The thermal shift assay is a rapid and inexpensive technique that quantifies change in protein denaturation temperature, and thus can be used to screen conditions that affect protein thermal stability, such as protein mutations, ligand binding, and buffer formulations (like pH, salts, detergents, and other additives). These assays are rapid (typically about 30 minutes) and are performed on a quantitative PCR system. The thermal shift method is compatible with high throughput screening and requires much less protein than methods such as circular dichroism and differential scanning calorimetry.

To learn more about GloMelt™ and see how it compares to competitor products, visit our GloMelt™ Technology Page.

Possible Uses for Thermal Shift Assays

- Optimize buffer formulation for protein stability and storage
- Determine how mutations affect your protein's stability
- Rapidly screen small molecule drug candidates and other ligands for protein binding

GloMelt™ Dye vs. Competitors

GloMelt™ dye has significant advantages over other environmentally sensitive dyes, such as SYPRO® Orange and PROTEOSTAT® TS dye. GloMeltTM dye generates a strong signal because it is optimized for detection in the SYBR® Green channel of qPCR instruments, and therefore low reaction volumes and low protein concentrations can be used. GloMelt™ dye is compatible with high concentrations of protein stabilizers (such as glycerol and sorbitol), and also protein destabilizers (such as DTT and imidazole). GloMelt™ dye performs very well in high detergent concentrations, unlike SYPRO® Orange. Another advantage is that ROX dye can be included with GloMelt™ dye during thermal shift assays, which improves results by increasing replicate consistency in PCR instruments that require ROX passive

SYBR is a registered trademark of Thermo Fisher Scientific. Proteostat is a registered trademark of Enzo Life Sciences.

This datasheet was generated on November 2, 2025 at 04:51:41 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/glomelt-thermal-shift-protein-stability-kit/