

**JBS Thermofluor Dye**

Cat. No.	Amount
X-TD-70	70 µl

**For general laboratory use.**

**Shipping:** shipped at ambient temperature

**Storage Conditions:** store at -20 °C

**Additional Storage Conditions:** protect from light

**Shelf Life:** 12 months

**Form:** liquid

**Concentration:** 5000x in DMSO

**Applications:**

JBS Thermofluor Dye complements JBScreen Thermofluor SPECIFIC HTS (CSS-333).

Fluorescence readout is recommended in the range of 570-600 nm.

**Related Products:**

JBScreen Thermofluor SPECIFIC HTS, #CS-333