

iFluor™ Labeling Products

Limited Time Only

**25%
OFF**

Use promo code: ifluor25

Coupon code ifluor25 must be provided at the time of sales for promotion to apply. Offer only applies to purchases of iFluor™ labeling products. Unless otherwise stated, offer only applies to purchases directly from AAT Bioquest, Inc. Contact us for purchases outside of USA. Offer expires on 5/31/2019.