



NEN® Life Science Products, Inc.  
Boston, MA 02118

## Technical Data Certificate of Analysis

Caution: For Laboratory Use. A research chemical for research purposes only.

### TSA™-Plus Fluorescence Palette System Catalog # NEL760

The TSA Plus Fluorescence Palette System contains four different fluorescently labeled tyramides and associated reagents; TSA Plus Fluorecein, TSA Plus TMR, TSA Plus Cyanine 3 and TSA Plus Cyanine 5. The kit contains enough reagents to stain 10-30 slides for evaluation of TSA technology.

**\*NOTE:** The enclosed manual should be followed EXCEPT when reconstituting the solid fluorophore tyramide in DMSO. For this procedure, please follow the directions below.

#### Stock Solution

##### **Fluorophore Tyramide (Amplification Reagent)**

**Fluorophore Tyramides (Amplification Reagents)** are supplied as a solid. Each vial must be reconstituted with 60 µL DMSO (dimethyl sulfoxide -molecular biology or HPLC-grade) before use. The Fluorophore Tyramide Stock Solution, when stored at 4° C, is stable for at least 3 months. (Note: DMSO freezes at 4° C. Thaw the Stock Solution before each use.)

#### Working Solution

##### **Fluorophore Tyramide(Amplification Reagent)**

Before each procedure, dilute the Fluorophore Tyramide Stock Solution 1:50 using **1X Plus Amplification Diluent** to make the Fluorophore Tyramide Working Solution. Approximately 100-300 µL of Fluorophore Tyramide Working Solution is required per slide. Discard any unused portion of working solution.