

## **Amido Black**

**Description:** **Amido** black is a versatile, inexpensive, protein dye for gels, PVDF, and nitrocellulose.

### **Protocols:**

#### **I Staining of acrylamide gels:**

Gels may be immersed in **amido** black working solution (1% w/v **amido** black in 7% acetic acid) for one hour and destained in changes of warm 3% acetic acid until clear.

#### **II Staining of Nitrocellulose Membranes<sup>1</sup>:**

Membranes are stained for 5 min in 40% **methanol/10%** acetic acid plus 0.5% (w/v) **amido** black, and then destained for 2 min in 40% methanol/10% in water.

**Note:** 0.01% **amido** black can also be **used<sup>2</sup>**.

#### **III Staining of PVDF:**

Membrane is stained in **amido** black working solution (0.1% **amido** black in 10% methanol, 2% acetic acid) for 10 to 15 min. Membranes are destained in 45% methanol, 7% acetic for 10 to 15 min.

To remove the last traces of stain, the membranes are immersed in 90% methanol and 7% acetic acid for 1 to 2 **mins**.

1. Harper et. al., ***Anal. Biochem.***, **157**, 270-274 (1986)

2. Sportsman, **J.R.** and Elder, J.H., ***Anal. Biochem.***, **139**, 298-300 (1984)