

NB100-129 Protocol

Western Blot Protocol for BMAL1 Antibody (NB100-129)

Procedure Guide for NB 100-129
(Anti-MOP3 aka Anti-BMAL1)

Western Blot Procedure

1. Proteins are separated by electrophoresis.
2. Proteins are transferred to nitrocellulose sheets.
3. Following the protein transfer, the filters are blocked for 1 hour at room temperature, gently shaking.
4. The filters are then incubated with a 1:100 dilution of (NB100-129) anti-MOP3 (aka: Banti-BMAL1) in PBS + 5% Dry Milk for 2 hours at room temperature, gently shaking. *
5. The filters are washed three times with TBST at room temperature.
6. The filters are then incubated with an Alkaline Phosphatase conjugated secondary antibody [Goat anti-Rabbit Alk Phos conjugated Ab from Jackson ImmunoResearch #111-055-003], diluted 1:5000 in PBS + 5% Dry Milk, for 1 hour at room temperature.
7. The filters are washed three times with TBST at room temperature.
8. Stained the filters with NBT/BCIP substrate at room temperature.
9. The filters are finally washed three times with TBST at room temperature.

* Overnight at 4 degrees C will increase sensitivity