



Orders: [orders@novusbio.com](mailto:orders@novusbio.com)

Support: [technical@novusbio.com](mailto:technical@novusbio.com)

Web: [www.novusbio.com](http://www.novusbio.com)

Protocols, Publications, Related Products, Reviews and more:

[www.novusbio.com/NB100-355](http://www.novusbio.com/NB100-355)

## NB100-355 Protocol

# Western Blot Protocol for RPE65 Antibody (NB100-355)

---

### Western Blot

1. Bovine RPE membrane (5ug) and recombinant human RPE65 (~20ug ) expressed in COS7 cells were loaded on a NuPAGE Novex Bis-Tris 10%gel along with SeeBlue Plus2 marker.
2. Run in NuPAGE MOPS electrophoresis buffer (Invitrogen).
- 3 Samples were probed with NB 100-355 RPE65 antibody (1:5,000 ) followed by alkaline phosphatase-conjugated goat anti-mouse IgG incubation.
4. Visualized by NBT/BCIP substrate.