



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/NB500-608](http://www.novusbio.com/NB500-608)

## NB500-608 Protocol

# Western Blot protocol for CD34 Antibody (NB500-608)

---

Western Blot Protocol for CD34 Antibody (NB500-608): [https://www.novusbio.com/products/cd34-antibody-4h11-apg\\_nb500-608](https://www.novusbio.com/products/cd34-antibody-4h11-apg_nb500-608)

Use Kg-1a human leukemia cell lysate as a positive control and JURKAT human leukemia T-cell line as a negative control.

Sample preparation: Resuspend approx. 50 million cells in 1 ml cold Lysis buffer (1% laurylmaltoside in 20 mM Tris/Cl, 100 mM NaCl pH 8.2, 50 mM NaF including Protease inhibitor Cocktail). Incubate 60 min on ice. Centrifuge to remove cell debris. Mix lysate with non-reducing SDS-PAGE sample buffer.

Application note: non-reducing conditions.