



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/NB100-123](http://www.novusbio.com/NB100-123)

## NB100-123 Protocol

### Immunohistochemistry protocol for HIF-1 alpha Antibody (NB100-123)

---

[[URL:[https://www.novusbio.com/products/hif-1-alpha-antibody-h1alpha67\\_nb100-123](https://www.novusbio.com/products/hif-1-alpha-antibody-h1alpha67_nb100-123)]][[Caption:HIF-1 alpha Antibody]]

Please see:

Primary Reference: Zhong, H., et al. Overexpression of Hypoxia-inducible Factor 1alpha in Common Human Cancers and Their Metastases. *Cancer Research*. 59: 5830-5835, 1999.