



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/NBC1-18526](http://www.novusbio.com/NBC1-18526)

## NBC1-18526 Protocol

### SDS-Page protocol for Serpin E1/PAI-1 Recombinant Protein (NBC1-18526)

SDS-PAGE (NBC1-18526): [https://www.novusbio.com/products/serpin-e1-pai-1-recombinant-protein\\_nbc1-18526](https://www.novusbio.com/products/serpin-e1-pai-1-recombinant-protein_nbc1-18526)

1. Prepare a recombinant PAI-1 with various concentrations (2,000 nM, 1,000 nM, 500 nM, 250 nM, 125 nM) in assay buffer. (Assay buffer: 50 mM Tris-HCl pH 8.0, 0.01% Tween 20)
2. Mix equal volumes of the PAI-1 protein and diluted 2 ug/ml uPA
3. Incubate at room temperature for 15 minutes.
4. After dilute incubated reaction mixtures (PAI-1, uPA) by a 5 fold dilution in assay buffer, mix 50ul of 200uM substrate (Z-GGR-AMC) and diluted mixtures.
4. Read excitation wavelength 355nm and emission wavelength 460nm for 10 minutes.