

PAD Cocktail, Active

Recombinant protein expressed in Sf9 cells

Catalog # P312-37C

Lot # Q2535-2

Product Description

The five full length recombined PAD proteins (1, 2, 3, 4 and 6) expressed by baculovirus in Sf9 insect cells using N-terminal GST tags. This mixture is ideally suited for in vitro citrullination of protein targets.

Molecular Weight

PAD1: approx. 95 kDa
PAD2: approx. 99 kDa
PAD3: approx. 97 kDa
PAD4: approx. 96 kDa
PAD6: approx. 100 kDa

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

Storage and Stability

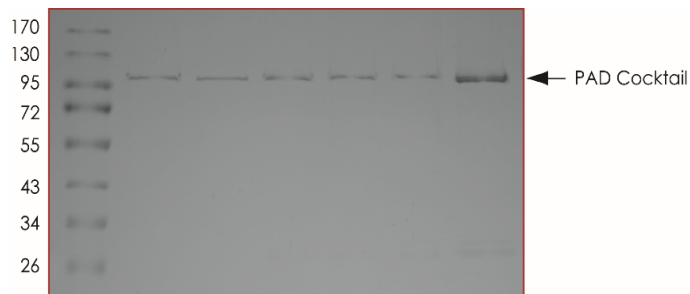
Store product at -70°C. For optimal storage, aliquot cocktail into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

Specific Activity

Sample Mass Spectrometry Data

Peptide Sequence	PAD Treated	Negative Control
XXRXXXXXXXXXX	5 reads, 3 citrullination	No citrullination
XXXXXXXXXXRXXRXXXXX	8 reads, 4 citrullination	No citrullination
XXXXXXXXRXXXXXXXXX	16 reads, 6 citrullination	No citrullination

Purity



SDS-PAGE image of PAD Cocktail Proteins. The five PAD isoform proteins from left to right are: PAD1, PAD4, PAD3, PAD6 and PAD2. The far right lane contains PAD Cocktail. The purity of PAD cocktail was determined to be **>95%** by densitometry, approx. MW **95-100 kDa**.

PAD Cocktail, Active

Recombinant protein expressed in Sf9 cells

Catalog #	P312-37C
Lot #	Q2535-2
Purity	>95%
Concentration	0.5 µg/µl (Each PAD at 0.1 µg/µl)
Stability	1yr at -70°C from date of shipment
Storage & Shipping	Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.

In Vitro Citrullination Protocol

Reaction Components

PAD Cocktail, Active (Catalog #: P312-37C)

10µl active PAD Cocktail (5µg of total PAD cocktail proteins) diluted with 20µl of PAD Cocktail Dilution Buffer (0.1M Tris-HCl, pH 7.4, 10mM CaCl₂) (Note: these are suggested working dilutions and it is recommended that the researcher perform appropriate dilutions for optimal results).

Anti-GST Agarose Beads

Any commercially available or self-made Anti-GST Agarose Beads may be used.

PAD Cocktail Dilution Buffer

Buffer components: 0.1M Tris-HCl, pH 7.4, 0.25mM DTT and 10mM CaCl₂.

PAD Buffer

Buffer components: 0.1M Tris-HCl, pH 7.4, 10mM CaCl₂ with 5mM DTT, 10µg/µl aprotinin, 10µg/µl leupeptin and 10µg/µl pepstatin.

4x Laemelli Sample Buffer

8% SDS, 40% glycerol, 250mM Tris-HCl (pH 6.8), 0.02% Bromophenol blue, and 8% β-mercaptoethanol.

Substrate Protein

Protein target to be determined by the researcher.

Citrullination Protocol

- Step 1.** Wash and equilibrate 5µl of anti-GST agarose beads in PAD Cocktail Dilution Buffer. Remove as much buffer as possible after equilibration.
- Step 2.** Incubate the equilibrated anti-GST agarose beads with 5µg of PAD Cocktail suspended in 20µl of PAD Cocktail Dilution Buffer at 4°C for 1 hour while shaking.
 - Component 1.** 5µl of equilibrated anti-GST agarose beads
 - Component 2.** 10µl of PAD Cocktail, Active (Catalog #P312-37C)
 - Component 3.** 20µl PAD Cocktail Dilution Buffer
- Step 3.** For a negative control, incubate the equilibrated anti-GST agarose beads in 20µl of PAD Cocktail Dilution Buffer.
- Step 4.** After incubation, wash the agarose beads three times in PAD buffer.
- Step 5.** Add 10µg of a Substrate Protein and 100µl PAD buffer to the PAD cocktail conjugated agarose beads and negative control, and incubate at 37°C for 2 hours while shaking.
 - Component 1.** 10µg of Substrate Protein
 - Component 2.** 100µl of PAD Buffer
- Step 6.** Centrifuge the beads and collect the supernatant containing the citrullinated Substrate Protein.
- Step 7.** Stop the reaction with 25µl of 4x Laemelli Sample Buffer.
- Step 8.** Store samples at -80°C until further use.
- Step 9.** To assess the level of citrullination and characterize the modification, the citrullinated and control samples can be sent for mass spectrometry analysis.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.