



Angiotensin II inhibition cocktail

Cat No: D05006 - for 10 ml of blood

General Data

Shipping:	dry ice
Stability:	Store at -80 degrees; shelf life 1 year maximum after production
Formulation:	Mix content of vial B into vial A (maximum volume of vial A is 10 mL) to obtain the ready-to-use 33x inhibition cocktail solution. Flocculation may appear.



Application(s): Prevent Angiotensin II formation and degradation

Product Overview

The Angiotensin II (All) Inhibition Cocktail has been developed by Bertin Bioreagent in order to fix physiological concentration of Angiotensin II in biological samples namely blood. Indeed, it has been observed that sample preparation is highly recommended when dosing Angiotensin II for preventing the formation or degradation of All in vitro.

Mix these solutions vial A and vial B to obtain the inhibitor cocktail ready to use. This solution has to be diluted 1:33 directly into the blood sample (i.e. for 1 mL of sample, add 30 μ L of inhibition cocktail). The total volume is for approximately one sample of 10 mL of blood. After adding the inhibition cocktail, mix the tube and centrifuge at 3000g for 15 minutes at +4°C.

Important notes: Plasma has to be prepared less than 30 minutes after collecting blood. Once mixed, the inhibition cocktail should be used immediately.

FP/27/19

For research laboratory use only - Not for human diagnostic use.

Buyers agree to purchase the material subject to Purchasing Terms that can be found on our website. Seek appropriate training to safely handle this product under normal conditions. Use the recommended personal protective equipment to prevent chemical exposures.

Bertin Bioreagent does not make any other warranty or representation whatsoever whether expressed or implied, with respect to these products. In no event will Bertin Bioreagent be liable for incidental, consequential or punitive damages.

Contact Bertin Bioreagent

Parc d'Activités du Pas du Lac
10 bis avenue Ampère
78180 Montigny le Bretonneux -
France

Tel.: +33 (0)139 306 036

<http://www.bertin-bioreagent.com>



Angiotensin II inhibition cocktail

Cat No: D05006 - for 50 ml of blood

General Data

Shipping:	dry ice
Stability:	Store at -80 degrees; shelf life 1 year maximum after production
Formulation:	Mix content of vial B into vial A (maximum volume of vial A is 10 mL) to obtain the ready-to-use 33x inhibition cocktail solution. Flocculation may appear.



Application(s): Prevent Angiotensin II formation and degradation

Product Overview

The angiotensin II (All) Inhibition Cocktail has been developed by Bertin Bioreagent in order to fix physiological concentration of Angiotensin II in biological samples namely blood. Indeed, it has been observed that sample preparation is highly recommended when dosing Angiotensin II for preventing the formation or degradation of All in vitro.

Mix these solutions vial A and vial B to obtain the inhibitor cocktail ready to use. This solution has to be diluted 1:33 directly into the blood sample (i.e. for 1 mL of sample, add 30 μ L of inhibition cocktail). The total volume is for approximately one sample of 50 mL of blood. After adding the inhibition cocktail, mix the tube and centrifuge at 3000g for 15 minutes at +4°C.

Important notes: Plasma has to be prepared less than 30 minutes after collecting blood. Once mixed, the inhibition cocktail should be used immediately.

FP/27/19

For research laboratory use only - Not for human diagnostic use.

Buyers agree to purchase the material subject to Purchasing Terms that can be found on our website. Seek appropriate training to safely handle this product under normal conditions. Use the recommended personal protective equipment to prevent chemical exposures.

Bertin Bioreagent does not make any other warranty or representation whatsoever whether expressed or implied, with respect to these products. In no event will Bertin Bioreagent be liable for incidental, consequential or punitive damages.

Contact Bertin Bioreagent

Parc d'Activités du Pas du Lac
10 bis avenue Ampère
78180 Montigny le Bretonneux -
France

Tel.: +33 (0)139 306 036

<http://www.bertin-bioreagent.com>



Angiotensin II inhibition cocktail

Cat No: D05006 - for 200 ml of blood

General Data

Shipping:	dry ice
Stability:	Store at -80 degrees; shelf life 1 year maximum after production
Formulation:	Mix content of vial B into vial A (maximum volume of vial A is 10 mL) to obtain the ready-to-use 33x inhibition cocktail solution. Flocculation may appear.



Application(s): Prevent Angiotensin II formation and degradation

Product Overview

The Angiotensin II (All) Inhibition Cocktail has been developed by Bertin Bioreagent in order to fix physiological concentration of Angiotensin II in biological samples namely blood. Indeed, it has been observed that sample preparation is highly recommended when dosing Angiotensin II for preventing the formation or degradation of All in vitro.

Mix these solutions vial A and vial B to obtain the inhibitor cocktail ready to use. This solution has to be diluted 1:33 directly into the blood sample (i.e. for 1 mL of sample, add 30 μ L of inhibition cocktail). The total volume is for approximately one sample of 200 mL of blood. After adding the inhibition cocktail, mix the tube and centrifuge at 3000g for 15 minutes at +4°C.

Important notes: Plasma has to be prepared less than 30 minutes after collecting blood. Once mixed, the inhibition cocktail should be used immediately.

FP/27/19

For research laboratory use only - Not for human diagnostic use.

Buyers agree to purchase the material subject to Purchasing Terms that can be found on our website. Seek appropriate training to safely handle this product under normal conditions. Use the recommended personal protective equipment to prevent chemical exposures.

Bertin Bioreagent does not make any other warranty or representation whatsoever whether expressed or implied, with respect to these products. In no event will Bertin Bioreagent be liable for incidental, consequential or punitive damages.

Contact Bertin Bioreagent

Parc d'Activités du Pas du Lac
10 bis avenue Ampère
78180 Montigny le Bretonneux -
France

Tel.: +33 (0)139 306 036

<http://www.bertin-bioreagent.com>