

# **OxyStat (BI-5007)**

## ***Frequently Asked Questions (FAQ)***

### ***Can the Oxystat kit be used on rat or mouse samples?***

The Oxystat kit is based on the reaction of peroxidase with plasma peroxides including lipid peroxides. So this is not specific for a certain species. Unfortunately we cannot provide you with a study using rat or mouse samples

### ***Which Plasma Peroxides are detected?***

The Oxystat kit is based on the reaction of peroxidase with plasma peroxides including lipid peroxides.  
So this is not specific for a certain species.