

生体イオウ関連研究用試薬

-SulfoBiotics- Series

