



BMG LABTECH

The Microplate Reader Company

Your research is too important to be left to others!



Come and see
our newest innovations
in microplate reading
technology at
LabAutomation '09,
booth 456

www.bmglabtech.com