

## Incident Log

Maxwell® RSC Instrument (Cat.# AS4500)

Record incidents of user or instrument error below.

Serial Number:

**Date of Incident** (MM/DD/YYYY): \_\_\_\_\_ **Time:** \_\_\_\_\_ **Ticket #:** \_\_\_\_\_

User Name: \_\_\_\_\_

Description of Incident: \_\_\_\_\_  
\_\_\_\_\_

Instrument Error Code (if shown): \_\_\_\_\_

Reported to Technical Services:  Yes /  No    If Yes, Scientist's Name: \_\_\_\_\_

Probable Cause: \_\_\_\_\_  
\_\_\_\_\_

Resolution: \_\_\_\_\_  
\_\_\_\_\_

Date Resolved (MM/DD/YYYY): \_\_\_\_\_

**Date of Incident** (MM/DD/YYYY): \_\_\_\_\_ **Time:** \_\_\_\_\_ **Ticket #:** \_\_\_\_\_

User Name: \_\_\_\_\_

Description of Incident: \_\_\_\_\_  
\_\_\_\_\_

Instrument Error Code (if shown): \_\_\_\_\_

Reported to Technical Services:  Yes /  No    If Yes, Scientist's Name: \_\_\_\_\_

Probable Cause: \_\_\_\_\_  
\_\_\_\_\_

Resolution: \_\_\_\_\_  
\_\_\_\_\_

Date Resolved (MM/DD/YYYY): \_\_\_\_\_