

dataTaker Power Transformer Change as of 26-03-2013

This document provides details of the change to a new switch mode power transformer for installation of <u>ALL</u> dataTaker products.

This Change applies to all users.

The transformer based power transformer is no longer available and has been replaced with an energy efficient switch mode power supply dataTaker Part Number PROMA-1000.

This global power supply has interchangeable input plug head, as shown below.

The power supply is MEPS and RoHS compliant, safety approvals UL/CUL/TUV/GS/SAA /CE/CB, safety compliant to UL 60950, EN 60950, AS NZS 60950 and IEC 60950.



IMPORTANT: When using the switch mode power supplies the dataTaker products <u>must</u> be Earthed / Grounded.

Please Note: General Safety Standards require that the logger is Earthed/Grounded.

An Earthing / Grounding point exists on all Loggers to meet these requirements.

Located on side panel as below – refer to Grounding section in product user manual.

