

Single Particle Optical Sizing

Preparation Checklist

- Suitable Working Area
- Location
- Electrical Services
- Rinse and Dilution Solution
- Drain Vessels
- Summary: Facilities Required



LPC 500 Liquid Particle Counter: *Standalone Particle Counting/ Sizing Configuration*

Introduction

The LPC 500™ liquid particle counter, Teledyne CETAC Oils 7400 autosampler and

the SPM 700 syringe module comprise a complete system for particle analysis of lubricating oils. A suitable working area, computer table or bench, and waste containers are still needed and must be provided by the analyst.

The complete particle counting system consists of the LPC 500 optical sensor, LPC 500 multi-channel pulse analyzer, Oils 7400 autosampler, SPM 700 syringe pump module, and the computer used to view the LPC 500 software interface to control the instrument – the dimensions of which are provided in Table 1 (Page 4).

