

## Solvent Compatibility and Utility

---

Most solvents typically used in life science laboratories are compatible with the fiber optic pedestals of all Thermo Scientific NanoDrop instruments. However, the high vapor pressure properties of some solvents may not be conducive to small volume measurements when using the pedestal for measurements on any of the NanoDrop™ Spectrophotometers. The cuvette option of the NanoDrop 2000/2000c Spectrophotometer is recommended when measuring samples with high vapor pressures.

The following solvents are compatible for use on the pedestals of all NanoDrop instruments: methanol, ethanol, n-propanol, isopropanol, butanol, acetone, ether, chloroform, carbon tetrachloride, DMSO, DMF, acetonitrile, THF, toluene, hexane, benzene, sodium hydroxide, sodium hypochlorite (bleach), dilute HCl, dilute HNO<sub>3</sub>, dilute acetic acid. It is recommended that all corrosive solvents be wiped from the pedestal immediately upon completion of a measurement. It is also recommended that the user end a series of measurements with a dH<sub>2</sub>O sample to ensure that solvents are not inadvertently left on the pedestal.

The diaphragm around the pedestal of the NanoDrop 2000/2000c is permanently affixed to the instrument. Do not attempt to remove the seal. Avoid prolonged exposure of the diaphragm to HCl, alcohol, bleach, acetone or other solvents as the adhesive securing the seal may be affected. If the seal comes loose please contact Technical Support.

All forms of Hydrofluoric Acid (HF) are incompatible as the fluoride ion will dissolve the quartz fiber optic cable.

For technical support, contact us at 302-479-7707 or [nanodrop@thermofisher.com](mailto:nanodrop@thermofisher.com).