

# Technical Bulletin

## TVA 2020 Cold Weather Operation

---

**Bulletin# TS0059****Rev12/15/15**

---

### Introduction

TVA2020 intrinsic safety design requires regulating the pump power to a level that can limit its ability to start at lower ambient temperatures (locked rotor condition). Attempting an FID ignition while the pump motor is in a locked rotor condition may trigger a safety circuit shutdown.

The TVA 2020 instruction manual specifies the “Normal Operating Conditions” as -10°C to +45°C. Due to this pump power limitation, the instrument low temperature start up condition is revised from -10°C to no less than 20°C. Once the pump is running, the operating temperatures are as originally specified (-10°C to +45°C).

Please be aware that a software resolution to this issue is under investigation and will be included in the next release, estimated completion in the first quarter of 2016.

### FID Ignition Procedure

- Option 1 - Store instrument in an environment and ignite above 20°C. Select 1=Run to start pump and ignite
- Option 2 - Start pump only (CONTROL 1) and let it run for 15 minutes prior to attempting to ignite