

# 3-A Antennas

## FOR VARIVENT® PIPELINE MOUNTING

### Robust & water resistant

IP65, designed for harsh operating conditions such as high temperatures, vibration, shock, and water. Maximum robustness against mechanical stress.

### Safety

Certified with 3-A Sanitary Standard

### Hygienic standard

CIP and SIP qualified.  
FDA suitable materials for the contact with food.



# TECHNICAL DATA & FACTS

## 3-A Antennas

### Antenna

---

Application	Food and beverages suitable for Varivent® pipe connectors
Temperature stability	Product temperature: 10 to 130 °C (283.to 403 K), temporarily up to 140 °C
Ambient temperature	-20 to +60 °C (253 to 333 K) operation -40 to +80 °C (283 to 353 K) storage
Connections	2 x HF connections: N female, 50 Ω for HF cable
Material	HF Antenna with a stainless-steel and PEEK microwave windows EPDM O-ring for fat/grease content lower than 8 % and HNBR O-ring if a fat/grease content higher than 8 %.
IP protection	IP 65
Certification	3-A Sanitary Standard(s) No.: 4066
Weight	1 kg

### Special Features

---

- Designed according to hygienic standards, CIP (Clean-In-Place) and SIP (Sterilize-In-Place) qualified, using FDA-compliant materials for food contact applications
- Certified with 3-A Sanitary Standard no. 4066
- Suitable for all Varivent® connectors