



## Signum® Complete Scale

Option Y2: for use in zone 2/22 hazardous areas (ATEX)



- For Signum Regular, Advanced, and Supreme scales
- For use in hazardous areas with potentially explosive gas and dust (non-conductive), in which an explosive atmosphere is not present, or is present only briefly, during normal operation (in acc. with Directive 1999/92/EU ATEX 137)
- Category 3GD equipment, designated and tested in accordance the new standard

The declaration of conformity for the Signum compact scale:



II 3 GD Ex nA nC nL IIC T4 T80°C IP6x

Manufacturer's declaration based on internal test report: SAG 06ATEX003X

- : Protection (dust tight, protected from penetration by jets of water)
- : All power supply options for the standard scale are available, as well as all accessories (except bar code scanner, Ethernet, below-scale weighing)
- : The manufacturer's declaration is made in accordance with the newest EN 60079-15 standard, edition 2005 (with regard to protection type "n"). The validity of the 2003 release with the previous "EEx" designation expires in 2007.
- : The built-in motorized calibration weight in the Signum Supreme (verified ex factory, standard equipment or optional) can also be used in conjunction with Option Y2
- : Permissible ambient operating temperature range: -10°C to +40°C (+14°F to +104°F)

