

AL-230

Technical Bulletin

Pressure Conversion Chart

To convert from	To	Multiply by	To convert from	To	Multiply by
bar	psi	14.5	micron (μ)	psi	0.00019337
bar	in. of H ₂ O	401.47	micron (μ)	in. of H ₂ O	0.0053525
bar	in. of Hg	29.53	micron (μ)	in. of Hg	0.0003937
bar	kPa	100	micron (μ)	kPa	0.0013332
bar	mbar	1000	micron (μ)	mbar	0.013332
bar	cm of H ₂ O	1019.73	micron (μ)	cm of H ₂ O	0.013595
bar	mm of Hg	750.06	micron (μ)	mm Hg	0.001
in. of H ₂ O	psi	0.036127	mm of Hg	psi	0.019337
in. of H ₂ O	in. of Hg	0.073554	mm of Hg	in. of H ₂ O	0.53525
in. of H ₂ O	kPa	0.2491	mm of Hg	in. of Hg	0.03937
in. of H ₂ O	mbar	2.491	mm of Hg	kPa	0.13332
in. of H ₂ O	cm of H ₂ O	2.54	mm of Hg	mbar	1.3332
in. of H ₂ O	mm of Hg	1.8683	mm of Hg	cm of H ₂ O	1.3595
in. of Hg	psi	0.4912	psi	in. of H ₂ O	27.68
in. of Hg	in. of H ₂ O	13.596	psi	in. of Hg	2.036
in. of Hg	kPa	3.3864	psi	kPa	6.8947
in. of Hg	mbar	33.864	psi	mbar	68.947
in. of Hg	cm of H ₂ O	34.532	psi	cm of H ₂ O	70.308
in. of Hg	mm of Hg	25.4	psi	mm of Hg	51.715
kPa	psi	0.14504	torr	psi	0.019337
kPa	in. of H ₂ O	4.0147	torr	in. of H ₂ O	0.53525
kPa	in. of Hg	0.2953	torr	in. of Hg	0.03937
kPa	mbar	10	torr	kPa	0.13332
kPa	cm of H ₂ O	10.1973	torr	mbar	1.3332
kPa	mm of Hg	7.5006	torr	cm of H ₂ O	1.3595
			torr	mm Hg	1
mbar	psi	0.0145			
mbar	in. of H ₂ O	0.40147			
mbar	in. of Hg	0.02953			
mbar	kPa	0.1			
mbar	cm of H ₂ O	1.01973			
mbar	mm of Hg	0.75006			