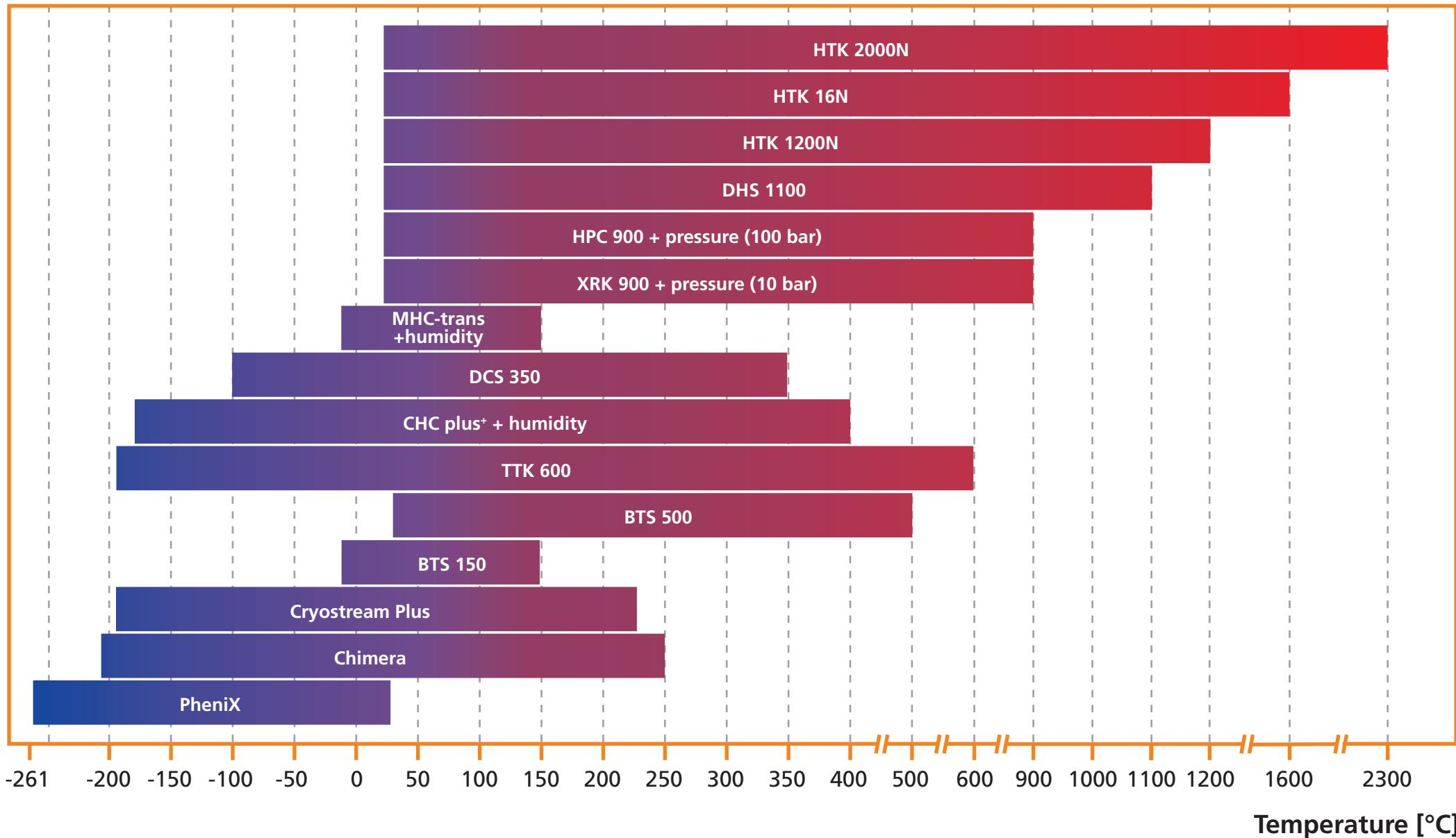


## Overview of non-ambient attachments (1)



# Overview of non-ambient attachments (2)

Attachment	Powder XRD	SAXS	PDF	GIXRD	Stress	XRR	Texture	High res. XRD	Air	Vacuum (mbar)	Inert gases	Humid gas	Reflection	Reflection + spin.	Transmission cap.	Reflection ChiPhi	Transmission foils
PheniX Cryostat	●			○	○	○				$10^{-4}$			●				
Chimera	●			○	○	○				$10^{-4}$			●				
Cryostream Plus	●	●	●						n/a	n/a	n/a	n/a			●		
BTS 150	○			○	○	○			●	$10^{-1}$	●		●				
BTS 500	○			○	○	○			●	$10^{-1}$	●		●				
TTK 600	●			○	○	○			●	$10^{-2}$	●		●		●		●
CHC plus+	●			○	○	○			●	$10^{-2}$	●	●	●				
DCS 350	○			●	●	●			●	$10^{-2}$	●					●	
MHC-trans	●	○							●		●	●					●
XRK 900	●			○	○	○			●		●		○	●			
HPC 900	●			○	○	○			●		●		●				
DHS 1100	○			●	●	●	●	●	●	$10^{-2}$	●					●	
HTK 1200N	●	○	○	○	○	○			●	$10^{-4}$	○		○	●	●		
HTK 16N	●								●	$< 10^{-4}$	●		●				
HTK 2000N	●								○	$< 10^{-4}$	●		●				

● = preferred    ○ = optional